

PRODUCT INFORMATION



Natural Products Screening Library (96-Well)

Item No. 20897 • Batch No. 0689479

Panels are routinely re-evaluated to include new catalog introductions as the research evolves.

Page 1 of 59

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Ursolic Acid	10072
1	A3	(+)-Absciscic Acid	10073
1	A4	(±)-Enterolactone	10112
1	A5	<i>trans</i> -trimethoxy Resveratrol	10188
1	A6	(R)-(-)-Mellein	10311
1	A7	Paclitaxel	10461
1	A8	Sulforaphane	10496
1	A9	Thapsigargin	10522
1	A10	Garcinol	10566
1	A11	Ellagic Acid	10569
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Apicidin	10575
1	B3	Isoliquiritigenin	10739
1	B4	(±)-Jasmonic Acid-Isoleucine	10740
1	B5	Epothilone B	10924
1	B6	Bisdemethoxycurcumin	10960
1	B7	Demethoxycurcumin	10961
1	B8	(R,S)-Anatabine	11001
1	B9	Piperlongumine	11006
1	B10	Delphinidin (chloride)	11012
1	B11	A23187	11016
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	all- <i>trans</i> Retinoic Acid	11017
1	C3	Forskolin	11018
1	C4	Bafilomycin A ₁	11038
1	C5	Betulin	11041
1	C6	Concanamycin A	11050
1	C7	4'-O-Methylhonokiol	11089
1	C8	Nicotinamide	11127
1	C9	Lupeol	11215
1	C10	Anisomycin	11308
1	C11	Ascomycin	11309
1	C12	<i>Unused</i>	

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/11/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 2 of 59

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Cyclopamine	11321
1	D3	Indole-3-carbinol	11325
1	D4	Cytochalasin A	11327
1	D5	Cytochalasin B	11328
1	D6	Cytochalasin D	11330
1	D7	Cytochalasin E	11331
1	D8	Fumagillin	11332
1	D9	Hypericin	11334
1	D10	Oligomycin A	11342
1	D11	Oligomycin B	11343
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Paxilline	11345
1	E3	Penitrem A	11347
1	E4	Withaferin A	11352
1	E5	Zearalenone	11353
1	E6	Mithramycin A	11434
1	E7	Mitomycin C	11435
1	E8	<i>Unused</i>	
1	E9	Sterigmatocystin	11441
1	E10	Tunicamycin Mixture	11445
1	E11	Phlorizin	11576
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	Natamycin	11634
1	F3	Amphotericin B	11636
1	F4	Pellitorine	11662
1	F5	12-Deoxywithastramonolide	11669
1	F6	14-deoxy-11,12-didehydro Andrographolide	11670
1	F7	α -acetyl Boswellic Acid	11673
1	F8	β -acetyl-Boswellic Acid	11674
1	F9	Aloe-emodin	11677
1	F10	Andrographolide	11679
1	F11	α -Asarone	11681
1	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	β -Asarone	11682
1	G3	Angelicin	11683
1	G4	Bakuchiol	11684
1	G5	Bavachin	11685
1	G6	Betulinic Acid	11686
1	G7	α -Boswellic Acid	11692
1	G8	Camptothecin	11694
1	G9	Catharanthine (tartrate)	11695
1	G10	Conessine	11700
1	G11	Deacetylforskolin	11701
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Demethylwedelolactone	11703
1	H3	6-Gingerol	11707
1	H4	Diacerein	11710
1	H5	β -Elemonic Acid	11712
1	H6	Gymnemagenin	11713
1	H7	Hexahydrocurcumin	11714
1	H8	Hypophyllanthin	11715
1	H9	Cycloastragenol	11722
1	H10	Jervine	11723
1	H11	Veratramine	11724
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	Platycodin D	11725
2	A3	Oleanolic Acid	11726
2	A4	Coumestrol	11730
2	A5	Ethyl <i>p</i> -methoxycinnamate	11740
2	A6	Neoandrographolide	11742
2	A7	Phyllanthin	11745
2	A8	Piperine	11750
2	A9	Purpurin	11752
2	A10	Rubiadin	11753
2	A11	Sesamolin	11754
2	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 4 of 59

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	<i>Unused</i>	
2	B3	(+)-ar-Turmerone	11759
2	B4	Vasicinone	11761
2	B5	Vinblastine (sulfate)	11762
2	B6	Vincamine	11763
2	B7	Vincristine (sulfate)	11764
2	B8	Vindoline	11765
2	B9	Vitamin D ₂	11791
2	B10	Vitamin D ₃	11792
2	B11	Wedelolactone	11796
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	Withanolide A	11797
2	C3	(-)-Epicatechin	11807
2	C4	(-)-Epicatechin Gallate	11808
2	C5	(-)-Epigallocatechin	11809
2	C6	4-Hydroxyisoleucine	11812
2	C7	Artemether	11815
2	C8	Artemisinin	11816
2	C9	Asiatic Acid	11818
2	C10	Asiaticoside	11819
2	C11	Bacopasaponin C	11820
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	Bacopaside I	11821
2	D3	Bacopaside II	11822
2	D4	Chebolic Acid	11829
2	D5	Colchicoside	11831
2	D6	Corilagin	11832
2	D7	Dioscin	11834
2	D8	Embelin	11838
2	D9	8-Gingerol	11841
2	D10	10-Gingerol	11842
2	D11	Glabridin	11843
2	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	18 β -Glycyrrhetic Acid	11845
2	E3	Gallic Acid	11846
2	E4	Glycyrrhizic Acid (ammonium salt)	11847
2	E5	Gracillin	11848
2	E6	Bacopaside X	11851
2	E7	Kaempferol	11852
2	E8	Licochalcone A	11853
2	E9	Madecassic Acid	11854
2	E10	Madecassoside	11855
2	E11	Brefeldin A	11861
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	Protodioscin	11887
2	F3	Rebaudioside A	11894
2	F4	Rebaudioside C	11895
2	F5	Atpenin A5	11898
2	F6	6-Shogaol	11901
2	F7	Stevioside (hydrate)	11902
2	F8	Trigonelline (chloride)	11904
2	F9	Salsalate	11911
2	F10	Ionomycin (calcium salt)	11932
2	F11	Triptolide	11973
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	Apocynin	11976
2	G3	Rottlerin	12006
2	G4	Atropine	12008
2	G5	Oxyresveratrol	12028
2	G6	Astragaloside IV	12069
2	G7	Etoposide	12092
2	G8	Kaempferitrin	12093
2	G9	Pterostilbene	13000
2	G10	Fucoxanthin	13068
2	G11	Punicalagin	13069
2	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 6 of 59

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	Pimecrolimus	13107
2	H3	Emodin	13109
2	H4	Anacardic Acid	13144
2	H5	(±)-Equol	13184
2	H6	cis-trismethoxy Resveratrol	13199
2	H7	Daidzin	13202
2	H8	trans-Zeatin	13226
2	H9	Rapamycin	13346
2	H10	Celastramycin A	13519
2	H11	Pseudolaric Acid B	13527
2	H12	<i>Unused</i>	
3	A1	<i>Unused</i>	
3	A2	Pristimerin	13621
3	A3	Icariin	13624
3	A4	Salidroside	13628
3	A5	Arecoline (hydrobromide)	13662
3	A6	Puromycin (hydrochloride)	13884
3	A7	Polydatin	13932
3	A8	Rotenone	13995
3	A9	Cafestol	13999
3	A10	Auraptene	14000
3	A11	(+)-Aphidicolin	14007
3	A12	<i>Unused</i>	
3	B1	<i>Unused</i>	
3	B2	Kahweol	14015
3	B3	Iberin	14016
3	B4	Erucin	14017
3	B5	Leupeptin (hemisulfate)	14026
3	B6	Ingenol	14031
3	B7	<i>Unused</i>	
3	B8	Caffeine	14118
3	B9	Indirubin	14155
3	B10	<i>Unused</i>	
3	B11	Glycitin	14161
3	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 7 of 59

Plate	Well	Contents	Item Number
3	C1	<i>Unused</i>	
3	C2	Glycitein	14162
3	C3	Naringenin	14173
3	C4	Genistin	14174
3	C5	Puerarin	14175
3	C6	α -Spinasterol	14197
3	C7	Magnolol	14233
3	C8	Tetrahydromagnolol	14234
3	C9	Febrifugine	14246
3	C10	Wogonin	14248
3	C11	Luotonin A	14253
3	C12	<i>Unused</i>	
3	D1	<i>Unused</i>	
3	D2	Luotonin F	14254
3	D3	Bilobalide	14272
3	D4	Hygromycin B (hydrate)	14291
3	D5	Plumbagin	14314
3	D6	Tanshinone IIA	14323
3	D7	Spectinomycin (hydrochloride hydrate)	14324
3	D8	<i>Unused</i>	
3	D9	Yangonin	14351
3	D10	Cordycepin	14426
3	D11	Asperlactone	14435
3	D12	<i>Unused</i>	
3	E1	<i>Unused</i>	
3	E2	Obscuroside A ₁	14439
3	E3	Tetrahydroharmine	14449
3	E4	Phloretin	14452
3	E5	Gossypol	14482
3	E6	<i>Unused</i>	
3	E7	Virginiamycin Complex	14503
3	E8	Ibotenic Acid	14584
3	E9	Cantharidin	14589
3	E10	Deferoxamine (mesylate)	14595
3	E11	Honokiol	14597
3	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 8 of 59

Plate	Well	Contents	Item Number
3	F1	<i>Unused</i>	
3	F2	7-dehydro Cholesterol	14612
3	F3	(-)-Huperzine A	14620
3	F4	Homoharringtonine	14631
3	F5	(S)-10-hydroxy-Camptothecin	14635
3	F6	Ginkgolide B	14636
3	F7	Aloesin	14649
3	F8	Xanthorrhizol	14668
3	F9	3-Nitropropionic Acid	14684
3	F10	Sophorose	14717
3	F11	1',4''-Sophorolactone 6',6''-diacetate	14718
3	F12	<i>Unused</i>	
3	G1	<i>Unused</i>	
3	G2	Costunolide	14743
3	G3	Cucurbitacin I	14747
3	G4	Shikonin	14751
3	G5	Gambogic Acid	14761
3	G6	Cyanidin (chloride)	14803
3	G7	Atractyloside (potassium salt)	14804
3	G8	Cucurbitacin B	14820
3	G9	Cucurbitacin E	14821
3	G10	Fusidic Acid (sodium salt)	14825
3	G11	Azaserine	14834
3	G12	<i>Unused</i>	
3	H1	<i>Unused</i>	
3	H2	<i>Unused</i>	
3	H3	<i>Unused</i>	
3	H4	Oxymatrine	14915
3	H5	Protocatechuic Acid	14916
3	H6	Digitonin	14952
3	H7	Polygodial	14979
3	H8	β -Lapachone	15021
3	H9	Thymoquinone	15039
3	H10	Phosphoramidon (sodium salt)	15113
3	H11	Vitexin	15116
3	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
4	A1	<i>Unused</i>	
4	A2	Psoralidin	15119
4	A3	(R,S)-Anatabine (tartrate)	15132
4	A4	Fisetin	15246
4	A5	A-54556A	15305
4	A6	(-)-Cotinine	15314
4	A7	Ginsenoside Rg ₁	15315
4	A8	Ginsenoside Rb ₁	15319
4	A9	<i>Unused</i>	
4	A10	Kasugamycin (hydrochloride)	15322
4	A11	Validamycin A	15326
4	A12	<i>Unused</i>	
4	B1	<i>Unused</i>	
4	B2	Ginsenoside Rd	15329
4	B3	Ginsenoside Re	15330
4	B4	Ginsenoside Rg ₂	15331
4	B5	20(S)-Ginsenoside Rg ₃	15332
4	B6	Harringtonine	15361
4	B7	Butein	15371
4	B8	Arctiin	15375
4	B9	Bicyclomycin Benzoate	15384
4	B10	Xanthohumol	15399
4	B11	Zerumbone	15400
4	B12	<i>Unused</i>	
4	C1	<i>Unused</i>	
4	C2	CAY10683	15403
4	C3	Nobiletin	15421
4	C4	Picroside I	15431
4	C5	Lasalocid	15505
4	C6	Phleomycin	15549
4	C7	Dobutamine (hydrochloride)	15582
4	C8	Elaiophylin	15583
4	C9	Rauwolscline (hydrochloride)	15596
4	C10	Daptomycin	15615
4	C11	Bufalin	15725
4	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
4	D1	<i>Unused</i>	
4	D2	Keracyanin (chloride)	15778
4	D3	3,3'-Diindolylmethane	15927
4	D4	Galangin	15948
4	D5	Vitamin K ₃	15950
4	D6	(R)-Prunasin	15959
4	D7	Fosmidomycin (sodium salt)	16000
4	D8	Roseoflavin	16001
4	D9	<i>Unused</i>	
4	D10	1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose	16007
4	D11	Tomatine	16024
4	D12	<i>Unused</i>	
4	E1	<i>Unused</i>	
4	E2	Muscone	16067
4	E3	Pheophorbide a	16072
4	E4	N-Acetylneuraminic Acid	16091
4	E5	Oenin	16092
4	E6	Delphinidin-3-β-D-glucoside (chloride)	16093
4	E7	DL-Sulforaphane N-acetyl-L-cysteine	16098
4	E8	<i>trans</i> -3'-Hydroxycotinine	16100
4	E9	Penta-N-acetylchitopentaose	16101
4	E10	<i>trans</i> -Clovamide	16138
4	E11	Sinomenine (hydrochloride)	16143
4	E12	<i>Unused</i>	
4	F1	<i>Unused</i>	
4	F2	Ingenol-3-angelate	16207
4	F3	Juglone	16216
4	F4	β-Nicotyrine	16333
4	F5	2,4-Diacetylphloroglucinol	16345
4	F6	Cyanidin 3-O-glucoside (chloride)	16406
4	F7	Reserpine	16474
4	F8	Biochanin A	16476
4	F9	Epothilone A	16482
4	F10	Erythromycin	16486
4	F11	Isorhamnetin	16496
4	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
4	G1	<i>Unused</i>	
4	G2	Spinosyn A	16528
4	G3	(±)-Nicotine	16535
4	G4	α-Zearalenol	16549
4	G5	Oleandomycin	16552
4	G6	Ferutinin	16554
4	G7	1-Methylnicotinamide (chloride)	16604
4	G8	3'-Sialyllactose (sodium salt)	16617
4	G9	Kanosamine (hydrochloride)	16623
4	G10	Monocrotaline	16666
4	G11	Astragaloside A	16677
4	G12	<i>Unused</i>	
4	H1	<i>Unused</i>	
4	H2	β-Carotene	16837
4	H3	8(Z),14(Z)-Eicosadienoic Acid	16854
4	H4	Swainsonine	16860
4	H5	(+)-Evodiamine	16885
4	H6	α-Naphthoflavone	16924
4	H7	Limonin	16932
4	H8	Tomatidine (hydrochloride)	16938
4	H9	(-)-Gallocatechin	16944
4	H10	7,8-Dihydroxyflavone	16946
4	H11	1,3,7-Trimethyluric Acid	16949
4	H12	<i>Unused</i>	
5	A1	<i>Unused</i>	
5	A2	Sanguinarine (chloride)	16951
5	A3	Indole-3-acetic Acid (sodium salt)	16954
5	A4	Casticin	16976
5	A5	Cryptotanshinone	16987
5	A6	A-54556B	17090
5	A7	Romidepsin	17130
5	A8	L-Sulforaphene	17152
5	A9	2,3,4,6-Tetra-O-benzoyl-β-D-glucopyranose	17163
5	A10	Sedanolid	17166
5	A11	<i>Unused</i>	
5	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
5	B1	<i>Unused</i>	
5	B2	Guvacoline (hydrobromide)	17190
5	B3	Noscapine (hydrochloride)	17255
5	B4	<i>Unused</i>	
5	B5	Rhein	17345
5	B6	Salicin	17357
5	B7	Phytol	17401
5	B8	Chrysin	17402
5	B9	Dihydrodaidzein	17423
5	B10	Evernic Acid	17430
5	B11	(5Z)-7-Oxozeaenol	17459
5	B12	<i>Unused</i>	
5	C1	<i>Unused</i>	
5	C2	8-Prenylnaringenin	17462
5	C3	Cytosporone B	17501
5	C4	Hinokitiol	17538
5	C5	Formononetin	17558
5	C6	Galantamine	17559
5	C7	Guanidinoethyl sulfonate	17572
5	C8	N-Acetyl-S-allyl-L-cysteine	17592
5	C9	<i>Unused</i>	
5	C10	N-(<i>p</i> -Coumaroyl) Serotonin	17797
5	C11	Salvianolic Acid A	17800
5	C12	<i>Unused</i>	
5	D1	<i>Unused</i>	
5	D2	DL-Cystathionine (hydrochloride)	17819
5	D3	6'-Sialyllactose (sodium salt)	17866
5	D4	Caffeic Acid methyl ester	17873
5	D5	Berteroin	17880
5	D6	N ⁶ -(Δ^2 -Isopentenyl)adenine	17906
5	D7	Myosmine	17908
5	D8	Tryptanthrin	17913
5	D9	Naringin	17923
5	D10	Micafungin (sodium salt)	18009
5	D11	Oxytetracycline	18076
5	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 13 of 59

Plate	Well	Contents	Item Number
5	E1	<i>Unused</i>	
5	E2	<i>Unused</i>	
5	E3	Capsiconiate	18172
5	E4	(±)-Anabasine (hydrochloride)	18206
5	E5	Digitoxigenin	18229
5	E6	Carbenoxolone (sodium salt)	18240
5	E7	<i>Unused</i>	
5	E8	Cardamonin	18310
5	E9	Anacardic Acid monoene	18422
5	E10	Menaquinone 4	18423
5	E11	Vitamin B ₁₂ (hydrate)	18425
5	E12	<i>Unused</i>	
5	F1	<i>Unused</i>	
5	F2	Galactinol (hydrate)	18426
5	F3	4-methyl-1-Pentanol	18446
5	F4	Boldine	18447
5	F5	all- <i>trans</i> Retinal	18449
5	F6	Patchouli Alcohol	18450
5	F7	Novobiocin (sodium salt)	18457
5	F8	Dehydrocostus lactone	18485
5	F9	Arphamenine B (hemisulfate)	18506
5	F10	Tyramine	18601
5	F11	Tropine	18627
5	F12	<i>Unused</i>	
5	G1	<i>Unused</i>	
5	G2	Gomisin C	18643
5	G3	(-)-Catechin	18644
5	G4	Quercetagetin	18645
5	G5	Hesperidin	18646
5	G6	(±)-Taxifolin	18647
5	G7	Quercetin 3-D-galactoside	18648
5	G8	Diosmetin	18649
5	G9	L-Histidinol (hydrochloride)	18739
5	G10	Ellipticine (hydrochloride)	18742
5	G11	Aureusimine B	18752
5	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
5	H1	<i>Unused</i>	
5	H2	Epothilone D	18765
5	H3	5-Phospho-D-ribose 1-diphosphate (sodium salt hydrate)	18897
5	H4	<i>Unused</i>	
5	H5	<i>Unused</i>	
5	H6	Mycophenolic Acid β -D-Glucuronide	19078
5	H7	Osthole	19195
5	H8	Syringic Acid	19196
5	H9	2-hydroxy-6-Methylbenzoic Acid	19199
5	H10	Thiostrepton	19200
5	H11	Nimbolide	19230
5	H12	<i>Unused</i>	
6	A1	<i>Unused</i>	
6	A2	<i>Unused</i>	
6	A3	Herbacetin	19285
6	A4	Esculetin	19286
6	A5	Bavachinin	19376
6	A6	Monoacetylphloroglucinol (hydrate)	19451
6	A7	Pleuromutilin	19452
6	A8	Clarithromycin	19455
6	A9	β -Zearalenol	19460
6	A10	Griseofulvin	19461
6	A11	Prodigiosin	19464
6	A12	<i>Unused</i>	
6	B1	<i>Unused</i>	
6	B2	<i>Unused</i>	
6	B3	Arglabin	19502
6	B4	<i>Unused</i>	
6	B5	Ferulenol	19568
6	B6	Podophyllotoxin	19575
6	B7	Sauchinone	19577
6	B8	1,7-Dimethyluric Acid	19584
6	B9	Eupatilin	19601
6	B10	Curvulin	19609
6	B11	Bafilomycin C ₁	19625
6	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
6	C1	<i>Unused</i>	
6	C2	6-Methylsulfinylhexyl isothiocyanate	19638
6	C3	β -Elemene	19641
6	C4	Juvenile Hormone III	19646
6	C5	Cepharanthine	19648
6	C6	D,L-1'-Acetoxychavicol Acetate	19649
6	C7	Racanisodamine	19650
6	C8	Dihydroresveratrol	19651
6	C9	Flavokawain B	19652
6	C10	<i>Unused</i>	
6	C11	Combretastatin A1	19750
6	C12	<i>Unused</i>	
6	D1	<i>Unused</i>	
6	D2	Sinensetin	19751
6	D3	Malvidin (chloride)	19752
6	D4	Pelargonidin (chloride)	19753
6	D5	Peonidin (chloride)	19754
6	D6	Petunidin (chloride)	19755
6	D7	7-methoxy Apigeninidin (chloride)	19757
6	D8	Jaceosidin	19758
6	D9	Vanillic Acid 4- β -D-glucopyranoside	19759
6	D10	6',7'-Epoxybergamottin	19760
6	D11	Sakuranetin	19761
6	D12	<i>Unused</i>	
6	E1	<i>Unused</i>	
6	E2	Z3-Dodecenyl E2-butenolate	19787
6	E3	Baicalin	19843
6	E4	Cinobufagin	19844
6	E5	Cinobufotalin	19845
6	E6	<i>Unused</i>	
6	E7	Diosmin	19848
6	E8	Schisandrin A	19849
6	E9	<i>Unused</i>	
6	E10	Isoimperatorin	19851
6	E11	Neobavaisoflavone	19852
6	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
6	F1	<i>Unused</i>	
6	F2	Mogroside V	19853
6	F3	Paeoniflorin	19861
6	F4	Paeonol	19862
6	F5	Physcion	19863
6	F6	Procyanidin B2	19865
6	F7	Quercitrin	19866
6	F8	Rhapontin	19867
6	F9	Rutin (hydrate)	19868
6	F10	Yohimbine (hydrochloride)	19869
6	F11	Chrysophanol	19870
6	F12	<i>Unused</i>	
6	G1	<i>Unused</i>	
6	G2	Ferulic Acid	19871
6	G3	Ginkgolide C	19872
6	G4	Isobavachalcone	19873
6	G5	Tetrandrine	19874
6	G6	7-hydroxy Coumarin Glucuronide (sodium salt)	19882
6	G7	Methyl Gallate	19951
6	G8	Loganin	19997
6	G9	6-Bromo-2-oxindole	20041
6	G10	Scopoletin	20042
6	G11	Norharmene	20043
6	G12	<i>Unused</i>	
6	H1	<i>Unused</i>	
6	H2	Trioxsalen	20044
6	H3	Homogentisic Acid	20045
6	H4	Scoparone	20046
6	H5	5-hydroxy-N ^ω -methyl Tryptamine (oxalate)	20047
6	H6	L-Quisqualic Acid	20211
6	H7	UCN-02	20221
6	H8	Icaritin	20236
6	H9	Vitamin A	20241
6	H10	Retinyl Acetate	20242
6	H11	Methysticin	20258
6	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
7	A1	<i>Unused</i>	
7	A2	16-O-Methylcafestol	20259
7	A3	Spiramycin	20267
7	A4	Rhamnetin	20302
7	A5	Epibrassinolide	20337
7	A6	Pyrogallol	20347
7	A7	<i>Unused</i>	
7	A8	<i>Unused</i>	
7	A9	Stachydrine (hydrochloride)	20506
7	A10	Folic Acid	20515
7	A11	Trehalose (hydrate)	20517
7	A12	<i>Unused</i>	
7	B1	<i>Unused</i>	
7	B2	N ⁶ -(Δ^2 -Isopentenyl)adenosine	20522
7	B3	Linamarin	20532
7	B4	L-Tetrahydropalmatine	20535
7	B5	Ansamitocin P-3	20538
7	B6	Sandramycin	20575
7	B7	Lumiflavine	20645
7	B8	Secologanin	20647
7	B9	5-(Hydroxymethyl)furfuryl Alcohol	20658
7	B10	Pyrroloquinoline quinone	20681
7	B11	Pyridoxine	20706
7	B12	<i>Unused</i>	
7	C1	<i>Unused</i>	
7	C2	7- <i>epi</i> Paclitaxel	20741
7	C3	N,N',N''-Triacetylchitotriose	20759
7	C4	Marinobufagenin	20798
7	C5	Isoquercitroside	20802
7	C6	Daphnetin	20826
7	C7	Acacetin	20827
7	C8	<i>Unused</i>	
7	C9	Nomilin	20836
7	C10	Homovanillic Acid sulfate (sodium salt)	20877
7	C11	(-)-Nicotine (tartrate)	20887
7	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
7	D1	<i>Unused</i>	
7	D2	Diallyl Sulfide	20894
7	D3	2-keto D-Glucose	20908
7	D4	Tyvelose	20910
7	D5	6-deoxy-L-Talose	20911
7	D6	Cephaeline	20916
7	D7	Brassinin	20922
7	D8	<i>p</i> -Coumaric Acid	20929
7	D9	3',4',7-Trihydroxyisoflavone	20930
7	D10	<i>Unused</i>	
7	D11	Tryptamine	20995
7	D12	<i>Unused</i>	
7	E1	<i>Unused</i>	
7	E2	β -Cembrenediol	21031
7	E3	Sisomicin	21046
7	E4	Emetine (hydrochloride hydrate)	21048
7	E5	Vitamin K ₁	21051
7	E6	3-Methylcytidine (methosulfate)	21064
7	E7	L-Iduronic Acid (sodium salt)	21066
7	E8	7-Methyluric Acid	21067
7	E9	Paraxanthine	21068
7	E10	Kukoamine B	21091
7	E11	Mesembrine	21100
7	E12	<i>Unused</i>	
7	F1	<i>Unused</i>	
7	F2	Carboxyatractyloside (potassium salt)	21120
7	F3	13-Methylberberine (chloride)	21154
7	F4	Isogarcinol	21164
7	F5	Bestatin	21217
7	F6	Oleuropein	21220
7	F7	Digallic Acid	21224
7	F8	3,4'-5-Trihydroxystilbene-3- β -D-Glucopyranoside	21246
7	F9	Quercetin 3-O-glucuronide	21289
7	F10	Pyropheophorbide <i>a</i>	21371
7	F11	<i>Unused</i>	
7	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
7	G1	<i>Unused</i>	
7	G2	Baccatin III	21404
7	G3	Tannic Acid	21421
7	G4	Glucotropaeolin (potassium salt)	21463
7	G5	1-Deoxynojirimycin	21500
7	G6	Aristolochic Acid	21520
7	G7	Lincomycin	21526
7	G8	Coenzyme B ₁₂ (hydrate)	21571
7	G9	β-Caryophyllene	21572
7	G10	(-)-Caryophyllene oxide	21573
7	G11	α-Humulene	21574
7	G12	<i>Unused</i>	
7	H1	<i>Unused</i>	
7	H2	(±)-Linalool	21575
7	H3	α-Pinene	21576
7	H4	(-)-β-Pinene	21577
7	H5	Dianemycin	21679
7	H6	Vulpinic Acid	21685
7	H7	9-cis Retinal	21692
7	H8	<i>Unused</i>	
7	H9	Terpinolene	21712
7	H10	Fascaplysin (chloride)	21715
7	H11	Mycophenolic Acid	21716
7	H12	<i>Unused</i>	
8	A1	<i>Unused</i>	
8	A2	(+)-Limonene	21719
8	A3	β-Myrcene	21720
8	A4	Theobromine	21745
8	A5	Curvularin	21758
8	A6	Demeclocycline (hydrochloride)	21759
8	A7	Amentoflavone	21779
8	A8	<i>Unused</i>	
8	A9	(-)-Perillyl Alcohol	21860
8	A10	Picrotin	21870
8	A11	Azadirachtin	21893
8	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
8	B1	<i>Unused</i>	
8	B2	(±)-Dunnione	21943
8	B3	Camalexin	21971
8	B4	Isochlortetracycline	22125
8	B5	(2'S)-Nicotine-1-oxide	22153
8	B6	Adipostatin A	22163
8	B7	Eucalyptol	22183
8	B8	Geranyl Acetate	22184
8	B9	<i>Unused</i>	
8	B10	D-Cycloserine	22194
8	B11	Timosaponin AIII	22236
8	B12	<i>Unused</i>	
8	C1	<i>Unused</i>	
8	C2	Aloperine	22259
8	C3	Cyclovirobuxine D	22260
8	C4	10-Deacetylbaicatin III	22261
8	C5	Cyanidin 3-O-β-D-Galactopyranoside (chloride)	22264
8	C6	Ginsenoside Rb ₂	22265
8	C7	Digoxin	22266
8	C8	Tubeimoside I	22276
8	C9	Elastatinal	22288
8	C10	Hydroquinidine	22295
8	C11	<i>Unused</i>	
8	C12	<i>Unused</i>	
8	D1	<i>Unused</i>	
8	D2	Mangiferin	22360
8	D3	Syringin	22387
8	D4	4'-O-methyl Quercetin	22406
8	D5	Procyanidin B1	22411
8	D6	Astragaloside I	22431
8	D7	Astragaloside III	22432
8	D8	Aloin	22435
8	D9	Pinostrobin	22436
8	D10	Brazilin	22456
8	D11	Biotin	22582
8	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 21 of 59

Plate	Well	Contents	Item Number
8	E1	<i>Unused</i>	
8	E2	Indole-3-butyric Acid	22591
8	E3	Urolithin A	22607
8	E4	Anacardic Acid diene	22663
8	E5	Orsellinic Acid	22708
8	E6	Ethyl Orsellinate	22709
8	E7	Kojic Acid	22712
8	E8	Isoferulic Acid	22715
8	E9	Incensole	22895
8	E10	Rutaecarpine	22897
8	E11	4'-hydroxy Chalcone	22898
8	E12	<i>Unused</i>	
8	F1	<i>Unused</i>	
8	F2	Matrine	22901
8	F3	Neohesperidin	23028
8	F4	Cyclophenin	23145
8	F5	Cardanol diene	23153
8	F6	Cardanol monoene	23154
8	F7	Cardanol triene	23155
8	F8	<i>p</i> -Cymene	23159
8	F9	(-)-Isopulegol	23161
8	F10	(+)-Fenchol	23162
8	F11	(R)-(+)-Pulegone	23163
8	F12	<i>Unused</i>	
8	G1	<i>Unused</i>	
8	G2	(+)-3-Carene	23165
8	G3	Geraniol	23166
8	G4	(+)-Valencene	23167
8	G5	(-)-Guaiol	23172
8	G6	Terpineol	23173
8	G7	(1R)-(+)-Camphor	23175
8	G8	(1S)-(-)-Camphor	23176
8	G9	Isoborneol	23177
8	G10	α -Terpinene	23178
8	G11	α -Phellandrene	23179
8	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
8	H1	<i>Unused</i>	
8	H2	Cedrene	23180
8	H3	Nerolidol	23182
8	H4	Sappanone A	23205
8	H5	Phloroglucinol	23223
8	H6	Acetosyringone	23224
8	H7	Isoxanthohumol	23233
8	H8	(+)-Taxifolin	23234
8	H9	Pseudomonic Acid	23238
8	H10	β -Methylamino-L-Alanine (hydrochloride)	23287
8	H11	Chlortetracycline	23302
8	H12	<i>Unused</i>	
9	A1	<i>Unused</i>	
9	A2	Aloin B	23321
9	A3	Guvacine (hydrochloride)	23361
9	A4	Cardol diene	23373
9	A5	Cardol triene	23374
9	A6	20(S)-Ginsenoside Rh ₂	23390
9	A7	Methylcardol triene	23404
9	A8	α -Zearalanol	23424
9	A9	Ginsenoside F ₂	23435
9	A10	(+)-Menthol	23463
9	A11	(+)- β -Citronellol	23464
9	A12	<i>Unused</i>	
9	B1	<i>Unused</i>	
9	B2	β -Ocimene	23466
9	B3	(+)-Borneol	23467
9	B4	(-)-Borneol	23468
9	B5	Deoxyviolacein	23477
9	B6	Brevianamide F	23493
9	B7	Terrein	23496
9	B8	Dihydromyricetin	23549
9	B9	Cellocidin	23637
9	B10	Ginsenoside Rf	23667
9	B11	γ -Terpinene	23722
9	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
9	C1	<i>Unused</i>	
9	C2	Retinyl Palmitate	23796
9	C3	Chloramphenicol Acetate	23811
9	C4	<i>Unused</i>	
9	C5	Phanerosporic Acid	23879
9	C6	Ergoloid (mesylates)	24095
9	C7	Hydroxocobalamin	24099
9	C8	Columbin	24105
9	C9	Transcroctetin	24217
9	C10	Diphenylamine	24250
9	C11	Nemorosone	24256
9	C12	<i>Unused</i>	
9	D1	<i>Unused</i>	
9	D2	(±)-Camphene	24478
9	D3	Phorbol	24479
9	D4	Umbelliferone	24665
9	D5	Voacangine	24692
9	D6	(-)-Securinine	24902
9	D7	Zingerone	24909
9	D8	(+)-Nootkatone	24910
9	D9	Atractylenolide III	24911
9	D10	Isosilybin	24913
9	D11	Amarogentin	24914
9	D12	<i>Unused</i>	
9	E1	<i>Unused</i>	
9	E2	(+)-Praeruptorin A	24915
9	E3	Gliorosein	24921
9	E4	Catalpol	24925
9	E5	Isoquercetin	24926
9	E6	Abietic Acid	24927
9	E7	Aucubin	24928
9	E8	Cyclo(D-Ala-L-Pro)	24940
9	E9	Cyclo(L-Leu-L-Pro)	24941
9	E10	Cyclo(L-Phe-L-Val)	24943
9	E11	Chicoric Acid	24960
9	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 24 of 59

Plate	Well	Contents	Item Number
9	F1	<i>Unused</i>	
9	F2	(-)-Gallocatechin Gallate	24962
9	F3	(-)-Catechin Gallate	24963
9	F4	3,5-Dicaffeoylquinic Acid	24964
9	F5	Verbascoside	24965
9	F6	Schizandrin B	24968
9	F7	Sennoside A	24969
9	F8	Ginkgolide A	24970
9	F9	Agnuside	24971
9	F10	<i>Unused</i>	
9	F11	Secoisolariciresinol Diglucoside	24974
9	F12	<i>Unused</i>	
9	G1	<i>Unused</i>	
9	G2	Maltotetraose	24975
9	G3	Notoginsenoside Ft ₁	24976
9	G4	Scutellarein	24977
9	G5	20(R)-Ginsenoside Rg ₃	24978
9	G6	4-deoxy-4 α -Phorbol	24998
9	G7	Maculosin	25019
9	G8	1,3-Dicaffeoylquinic Acid	25022
9	G9	Galactitol	25023
9	G10	Saikosaponin A	25026
9	G11	Aconine	25048
9	G12	<i>Unused</i>	
9	H1	<i>Unused</i>	
9	H2	Marmesin	25049
9	H3	1,5-Dicaffeoylquinic Acid	25050
9	H4	15,16-Dihydrotanshinone I	25051
9	H5	20(R)-Protopanaxadiol	25052
9	H6	3,4-Dicaffeoylquinic Acid	25053
9	H7	5-O-Caffeoylquinic Acid	25054
9	H8	Alkannin	25058
9	H9	Artemisinic Acid	25059
9	H10	Kaempferol-3-glucoside	25060
9	H11	Atractylenolide I	25061
9	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 25 of 59

Plate	Well	Contents	Item Number
10	A1	<i>Unused</i>	
10	A2	Baohuoside I	25062
10	A3	β -Mangostin	25071
10	A4	Caftaric Acid	25090
10	A5	Calycosin	25091
10	A6	Calycosin 7-O- β -D-Glucoside	25095
10	A7	Chelidonine	25097
10	A8	Cryptochlorogenic Acid	25098
10	A9	Cynaropicrin	25099
10	A10	Deacetylasperulosidic Acid	25100
10	A11	Dehydroevodiamine (hydrochloride)	25101
10	A12	<i>Unused</i>	
10	B1	<i>Unused</i>	
10	B2	Gentiopicrin	25102
10	B3	Ginkgetin	25103
10	B4	Isoginkgetin	25104
10	B5	(-)-Isoxanthohumol	25105
10	B6	3-methyl Orsellinic Acid	25113
10	B7	Cinnamamide	25117
10	B8	Magnoflorine	25122
10	B9	Magnolin	25123
10	B10	Narirutin	25124
10	B11	(+)-Gallocatechin	25125
10	B12	<i>Unused</i>	
10	C1	<i>Unused</i>	
10	C2	23- <i>epi</i> -26-Deoxyactein	25126
10	C3	Parishin C	25127
10	C4	Parishin E	25128
10	C5	Theaflavin	25129
10	C6	Ginsenoside Ro	25137
10	C7	Hederacoside C	25138
10	C8	Kukoamine A	25139
10	C9	Soyasaponin Bb	25140
10	C10	Luteoloside	25141
10	C11	Dihydroartemisinic Acid	25142
10	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
10	D1	<i>Unused</i>	
10	D2	Orientin	25143
10	D3	Kinsenoside	25144
10	D4	Senecionine	25145
10	D5	Seneciphylline	25146
10	D6	Ginsenoside Rg ₅	25147
10	D7	Lonicerin	25148
10	D8	Schisandrol B	25149
10	D9	Didymin	25197
10	D10	Isosakuranetin	25198
10	D11	Senkyunolide A	25199
10	D12	<i>Unused</i>	
10	E1	<i>Unused</i>	
10	E2	Procyanidin A2	25200
10	E3	Procyanidin C1	25201
10	E4	Chebulagic Acid	25203
10	E5	Voacamine	25204
10	E6	4-Hydroxy-6-methyl-2-pyrone	25210
10	E7	Decursinol angelate	25212
10	E8	Nodakenin	25213
10	E9	Decursin	25214
10	E10	Theaflavin 3,3'-digallate	25215
10	E11	Periplocin	25216
10	E12	<i>Unused</i>	
10	F1	<i>Unused</i>	
10	F2	Ruscogenin	25217
10	F3	Cryogenine	25287
10	F4	Thiamine (nitrate)	25332
10	F5	N-Acetylpuromycin	25335
10	F6	Micheliolide	25402
10	F7	(-)-Voacangarine	25468
10	F8	Dihydroaeruginic Acid	25514
10	F9	Mitraphylline	25528
10	F10	<i>Unused</i>	
10	F11	Thiamine (hydrochloride)	25656
10	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
10	G1	<i>Unused</i>	
10	G2	Oridonin	25665
10	G3	Benzoylaconine	25687
10	G4	Obafluorin	25731
10	G5	<i>trans</i> - ϵ -Viniferin	25737
10	G6	Piperitenone	25752
10	G7	($\pm$)-Menthol	25753
10	G8	($\pm$)-3-Carene	25768
10	G9	(+)-Cedrol	25769
10	G10	α -Copaene	25770
10	G11	β -Eudesmol	25771
10	G12	<i>Unused</i>	
10	H1	<i>Unused</i>	
10	H2	($\pm$)-Limonene	25773
10	H3	Nerol	25774
10	H4	Sabinene	25777
10	H5	<i>trans</i> -Nerolidol	25778
10	H6	Lythrine	25804
10	H7	Trilobatin	25837
10	H8	Neoeriocitrin	25839
10	H9	Lythridine	25865
10	H10	Hecogenin Acetate	25977
10	H11	Sinigrin (hydrate)	26028
10	H12	<i>Unused</i>	
11	A1	<i>Unused</i>	
11	A2	Dendrobine	26067
11	A3	Imperatorin	26070
11	A4	<i>Unused</i>	
11	A5	Corynecin V	26167
11	A6	Kirenol	26212
11	A7	(E)-Ajoene	26242
11	A8	Bergapten	26254
11	A9	Olivetolic Acid	26282
11	A10	Varinolic Acid	26283
11	A11	($\pm$)-Goitrin	26382
11	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
11	B1	<i>Unused</i>	
11	B2	Hispidulin	26383
11	B3	Specnuezhenide	26385
11	B4	Bergenin	26406
11	B5	Arbutin	26407
11	B6	Daphnoretin	26408
11	B7	Totarol	26412
11	B8	Peimisine (hydrochloride)	26438
11	B9	Lateropyrone	26463
11	B10	(+)-Kavain	26524
11	B11	<i>Unused</i>	
11	B12	<i>Unused</i>	
11	C1	<i>Unused</i>	
11	C2	Germacrene D	26539
11	C3	Agistatin B	26600
11	C4	Anacardic Acid triene	26611
11	C5	Parishin A	26665
11	C6	Amygdalin	26668
11	C7	Ganoderic Acid A	26674
11	C8	γ -Mangostin	26675
11	C9	Antipain (hydrochloride)	26705
11	C10	Curcumenol	26715
11	C11	Schisantherin B	26716
11	C12	<i>Unused</i>	
11	D1	<i>Unused</i>	
11	D2	($\pm$)- β -Tocopherol	26758
11	D3	Ganoderic Acid D	26765
11	D4	2S-Eriodictyol	26803
11	D5	Apigenin 7-glucoside	26813
11	D6	D-Pinitol	26815
11	D7	Vitexin 2-O-rhamnoside	26816
11	D8	Vatiquinone	26819
11	D9	Tabersonine	26848
11	D10	Curcumol	26849
11	D11	Leonurine	26850
11	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 29 of 59

Plate	Well	Contents	Item Number
11	E1	<i>Unused</i>	
11	E2	Shikimic Acid	26851
11	E3	Tanshinone I	26852
11	E4	Isoorientin	26862
11	E5	Farnesene (mixed isomers)	26891
11	E6	Linarin	26900
11	E7	Myricetrin	26902
11	E8	(Z)-Ligustilide	26908
11	E9	Altenusin	26911
11	E10	Ononin	27046
11	E11	<i>Unused</i>	
11	E12	<i>Unused</i>	
11	F1	<i>Unused</i>	
11	F2	(S)-Laudanosine	27170
11	F3	(±)-γ-Tocopherol	27193
11	F4	Schizandrin	27211
11	F5	Troxerutin	27220
11	F6	Brucine	27221
11	F7	Danshensu (sodium salt)	27222
11	F8	4β-Hydroxywithanolide E	27245
11	F9	Streptazolin	27266
11	F10	Ferrioxamine E	27275
11	F11	Complanatuside	27295
11	F12	<i>Unused</i>	
11	G1	<i>Unused</i>	
11	G2	Plantamajoside	27296
11	G3	20(R)-Ginsenoside Rh ₂	27300
11	G4	Jatrorrhizine (chloride)	27314
11	G5	Ginsenoside Rh ₁	27318
11	G6	Oroxylin A	27363
11	G7	D-Mannose	27388
11	G8	Nicotiflorin	27392
11	G9	Ligustroflavone	27431
11	G10	23-hydroxy Betulinic Acid	27446
11	G11	Loureirin B	27454
11	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
11	H1	<i>Unused</i>	
11	H2	Asarinin	27456
11	H3	N- <i>trans</i> -Feruloyltyramine	27459
11	H4	Santonin	27460
11	H5	Scutellarin	27461
11	H6	4-Hydroxybenzoic Acid	27468
11	H7	Schisandrin C	27475
11	H8	Echinocystic Acid	27476
11	H9	β -Rubromycin	27477
11	H10	Linderane	27479
11	H11	Aurantioobtusin	27481
11	H12	<i>Unused</i>	
12	A1	<i>Unused</i>	
12	A2	Hederoside D ₂	27507
12	A3	Mulberroside A	27545
12	A4	3,4-Dihydroxybenzaldehyde	27547
12	A5	6-Paradol	27554
12	A6	Dieugenol	27557
12	A7	Thymohydroquinone	27558
12	A8	Epimedin A	27582
12	A9	Monotropein	27595
12	A10	Acetylcimigenol Arabinoside	27596
12	A11	Tyrosol	27600
12	A12	<i>Unused</i>	
12	B1	<i>Unused</i>	
12	B2	Nepodin	27611
12	B3	Pinosylvin	27612
12	B4	Gentisic Acid	27615
12	B5	Questiomycin A	27623
12	B6	Oxychlororaphine	27624
12	B7	Swertiamarin	27634
12	B8	Quercetin-7-O- β -D-Glucopyranoside	27641
12	B9	Dihydromethysticin	27644
12	B10	Corydaline	27654
12	B11	Ginkgoneolic Acid	27656
12	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
12	C1	<i>Unused</i>	
12	C2	Carubicin	27660
12	C3	Columbianadin	27661
12	C4	Agarotetrol	27662
12	C5	Ganoderenic Acid A	27685
12	C6	(+)-Pinoresinol	27700
12	C7	Pinoresinol Diglucoside	27708
12	C8	Talatisamine	27721
12	C9	Narcissoside	27759
12	C10	Naringin Dihydrochalcone	27791
12	C11	Xanthotoxol	27801
12	C12	<i>Unused</i>	
12	D1	<i>Unused</i>	
12	D2	Phenethyl isothiocyanate	27804
12	D3	Betulinic Aldehyde	27810
12	D4	Digitoxin	27825
12	D5	Diphyllin	27826
12	D6	5-Hydroxymethylfurfural	27835
12	D7	Cephalomannine	27878
12	D8	Rebaudioside D	27880
12	D9	Corosolic Acid	27887
12	D10	4,5-Dicaffeoylquinic Acid	27892
12	D11	Umbelliprenin	27894
12	D12	<i>Unused</i>	
12	E1	<i>Unused</i>	
12	E2	Bilobetin	27899
12	E3	Sciadopitysin	27900
12	E4	Schaftoside	27950
12	E5	<i>Unused</i>	
12	E6	Cyclo(L-Pro-L-Val)	27961
12	E7	Astringin	27979
12	E8	Telocinobufagin	27981
12	E9	Rubimaillin	27999
12	E10	Fraxetin	28011
12	E11	1,6-Dimethoxyphenazine	28061
12	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 32 of 59

Plate	Well	Contents	Item Number
12	F1	<i>Unused</i>	
12	F2	Verrucofortune	28062
12	F3	2-O-Acetyl-20-hydroxyecdysone	28063
12	F4	1-Hydroxyphenazine	28064
12	F5	Atranorin	28106
12	F6	Solamargine	28110
12	F7	Solasonine	28111
12	F8	<i>Unused</i>	
12	F9	Garcinoic Acid	28122
12	F10	α -Hederin	28164
12	F11	Fraxin	28217
12	F12	<i>Unused</i>	
12	G1	<i>Unused</i>	
12	G2	Fraxinol	28218
12	G3	Crotonoside	28219
12	G4	Astilbin	28267
12	G5	Geniposide	28281
12	G6	Oxysophocarpine	28305
12	G7	($\pm$)- α -Tocopherol Acetate	28399
12	G8	Loganic Acid	28402
12	G9	Tryptophol	28404
12	G10	Geniposidic Acid	28409
12	G11	Sennoside B	28410
12	G12	<i>Unused</i>	
12	H1	<i>Unused</i>	
12	H2	Coptisine (chloride)	28424
12	H3	Nicotinamide N-oxide	28441
12	H4	Maltopentaose	28456
12	H5	Precocene II	28480
12	H6	Lupulone	28481
12	H7	Corynoxine (hydrochloride)	28496
12	H8	10-deacetyl Paclitaxel	28503
12	H9	Aristolochic Acid B	28504
12	H10	Aristolochic Acid C	28505
12	H11	Esculin (hydrate)	28514
12	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
13	A1	<i>Unused</i>	
13	A2	Hypocrellin B	28544
13	A3	Pulchrenoside A	28545
13	A4	Quercetin 3-O-sophoroside	28592
13	A5	Demethylzeylasteral	28595
13	A6	Paclitaxel octadecanedioate	28635
13	A7	SAHO	28725
13	A8	Lappaconitine	28747
13	A9	Aescin	28815
13	A10	Indole-3-propionic Acid	28821
13	A11	(±)-Liquiritigenin	28823
13	A12	<i>Unused</i>	
13	B1	<i>Unused</i>	
13	B2	Tenuigenin	28824
13	B3	Gastrodin	28828
13	B4	Panaxatriol	28835
13	B5	Rhein-8-glucoside	28844
13	B6	Morusin	28857
13	B7	Kaurenoic Acid	28957
13	B8	Ginsenoside Rb ₃	29005
13	B9	5,7-Dihydroxychromone	29006
13	B10	Rehmannioside D	29034
13	B11	Melittoside	29035
13	B12	<i>Unused</i>	
13	C1	<i>Unused</i>	
13	C2	Rubusoside	29040
13	C3	Forsythoside B	29051
13	C4	Procyanidin B3	29065
13	C5	20(S)-Protopanaxatriol	29073
13	C6	2''-O-Galloylhyperin	29086
13	C7	Ginsenoside Rc	29088
13	C8	4'-O-Methylresveratrol	29106
13	C9	α-Mangostin	29108
13	C10	Methylcobalamin	29113
13	C11	(-)-Nicotine	29138
13	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 34 of 59

Plate	Well	Contents	Item Number
13	D1	<i>Unused</i>	
13	D2	Cyclosporin C	29152
13	D3	(-)-Tryptonide	29156
13	D4	Ginkgolide J	29163
13	D5	(±)-Catechin (hydrate)	29175
13	D6	Polyphyllin I	29179
13	D7	2-heptyl-3-hydroxy-4(1H)-Quinolone	29186
13	D8	Ailanthone	29194
13	D9	Agistatin E	29211
13	D10	Isoverbascoside	29216
13	D11	Nitidine (chloride)	29223
13	D12	<i>Unused</i>	
13	E1	<i>Unused</i>	
13	E2	Norwogonin	29227
13	E3	Maytansinol	29228
13	E4	Fangchinoline	29243
13	E5	Gibberellic Acid	29322
13	E6	Diallyl Tetrasulfide	29328
13	E7	Wogonoside	29332
13	E8	Citropten	29338
13	E9	<i>trans</i> -β-Ocimene	29393
13	E10	Elemicin	29394
13	E11	Prunetin	29432
13	E12	<i>Unused</i>	
13	F1	<i>Unused</i>	
13	F2	Arjunolic Acid	29438
13	F3	Asperuloside	29439
13	F4	Pulsatilla Saponin D	29455
13	F5	Coniferyl Alcohol	29470
13	F6	(-)-Lobelin (hydrochloride)	29526
13	F7	Sophocarpine	29598
13	F8	Eupatorin	29599
13	F9	Josamycin	29606
13	F10	Harmane	29613
13	F11	(±)-Bornyl Acetate	29617
13	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
13	G1	<i>Unused</i>	
13	G2	Peiminine	29618
13	G3	Coniferyl Ferulate	29620
13	G4	Daurisoline	29628
13	G5	Isorhynchophylline	29650
13	G6	Phillyrin	29654
13	G7	Hypaconitine	29656
13	G8	Ligustrazine (hydrochloride)	29658
13	G9	Kaempferide	29673
13	G10	Tropolone	29674
13	G11	Eleutheroside E	29684
13	G12	<i>Unused</i>	
13	H1	<i>Unused</i>	
13	H2	Purpurogallin	29689
13	H3	Notoginsenoside R1	29691
13	H4	Epimedin B	29694
13	H5	13-acetyl-9-Dihydrobaccatin III	29701
13	H6	Gypenoside IX	29705
13	H7	Malabaricone C	29741
13	H8	Albiflorin	29744
13	H9	Alantolactone	29762
13	H10	Procyanidin	29763
13	H11	Oleanolic Acid Acetate	29764
13	H12	<i>Unused</i>	
14	A1	<i>Unused</i>	
14	A2	Betulonic Acid	29765
14	A3	Estragole	29824
14	A4	Pinocembrin	29852
14	A5	Homoplantagin	29853
14	A6	(E)-Anethol	29862
14	A7	Oleandrin	29871
14	A8	(S)-(-)-Perillaldehyde	29872
14	A9	Indole-3-pyruvic Acid	29876
14	A10	7,3',4'-Trihydroxyflavone	29880
14	A11	(-)-Isocorypalmine	29886
14	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
14	B1	<i>Unused</i>	
14	B2	(-)-Sinoacutine	29889
14	B3	Malabaricone B	29911
14	B4	Geraniin	29924
14	B5	Ginsenoside CK	29926
14	B6	Echinacoside	29927
14	B7	Hypocrellin A	29933
14	B8	Bruceine A	29936
14	B9	Ajmaline	29962
14	B10	Peimine	29963
14	B11	D-threo-Biopterin	29986
14	B12	<i>Unused</i>	
14	C1	<i>Unused</i>	
14	C2	5-Heneicosylresorcinol	30031
14	C3	5-Tricosylresorcinol	30032
14	C4	Deltorpin C (trifluoroacetate salt)	30050
14	C5	Australine (hydrochloride)	30051
14	C6	Polygalasaponin F	30064
14	C7	2-Hydroxy-4-methoxybenzaldehyde	30091
14	C8	Isocorydine	30101
14	C9	Myricetin 3,4'-di-O- β -glucopyranoside	30102
14	C10	Schisanhenol	30109
14	C11	Forsythoside A	30110
14	C12	<i>Unused</i>	
14	D1	<i>Unused</i>	
14	D2	Germacrone	30111
14	D3	Ganoderic Acid B	30112
14	D4	Neotuberostemonine	30116
14	D5	Dihydrochelerythrine	30118
14	D6	Erythrodiol	30135
14	D7	<i>Unused</i>	
14	D8	Lapachol	30190
14	D9	Methylophiopogonanone A	30192
14	D10	Methylophiopogonanone B	30193
14	D11	Gomisin J	30213
14	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
14	E1	<i>Unused</i>	
14	E2	Araloside A	30219
14	E3	Pseudoginsenoside F ₁₁	30222
14	E4	Pelargonidin 3-O-glucoside (chloride)	30224
14	E5	Liquiritin	30263
14	E6	Secoxyloganin	30267
14	E7	Isoliquiritin	30270
14	E8	Pulchrenoside B ₄	30271
14	E9	2,3,5,6-Tetramethylpyrazine	30281
14	E10	DL-3-n-Butylphthalide	30284
14	E11	Sclareol	30287
14	E12	<i>Unused</i>	
14	F1	<i>Unused</i>	
14	F2	Palmatine (chloride)	30318
14	F3	Picroside II	30319
14	F4	Pteryxin	30320
14	F5	Coniferin	30364
14	F6	Curdione	30365
14	F7	Ganoderenic Acid C	30366
14	F8	Ganoderic Acid C2	30371
14	F9	Lucidenic Acid A	30372
14	F10	(-)-Menthol	30384
14	F11	Tenuifolin	30617
14	F12	<i>Unused</i>	
14	G1	<i>Unused</i>	
14	G2	Polygalacic Acid	30620
14	G3	Polygalaxanthone III	30621
14	G4	Chrysosplenetin	30625
14	G5	Methyl 3,4-Dihydroxyphenylacetate	30637
14	G6	Hederasaponin B	30641
14	G7	Timosaponin BII	30644
14	G8	Lycorine	30650
14	G9	Montbretin A	30651
14	G10	Arenobufagin	30695
14	G11	Tormentic Acid	30696
14	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
14	H1	<i>Unused</i>	
14	H2	Gomisin D	30697
14	H3	Phillygenin	30698
14	H4	Quercetin 3-O- α -L-arabinopyranoside	30707
14	H5	Isocurcumenol	30710
14	H6	Morroniside	30711
14	H7	Corynoline	30712
14	H8	<i>trans</i> -Zeatin Riboside	30747
14	H9	Idronoxil	30783
14	H10	2(E),4(Z)-Decadienoic Acid ethyl ester	30835
14	H11	7-Methoxycoumarin	30857
14	H12	<i>Unused</i>	
15	A1	<i>Unused</i>	
15	A2	Ergosterol Peroxide	30860
15	A3	<i>Unused</i>	
15	A4	Deslanoside	30872
15	A5	Cirsiliol	30874
15	A6	Ethamivan	30876
15	A7	Brusatol	30883
15	A8	Fuziline	30889
15	A9	Eurycomanone	30896
15	A10	Mahanimbine	30897
15	A11	Denudatine	30904
15	A12	<i>Unused</i>	
15	B1	<i>Unused</i>	
15	B2	Garcinone D	30906
15	B3	(3R,8S)-Falcarindiol	30919
15	B4	Blinin	30945
15	B5	Dehydrocorydaline (chloride)	30972
15	B6	<i>Unused</i>	
15	B7	Afzelin	31135
15	B8	Ginsenoside F ₁	31137
15	B9	Dihydrosanguinarine	31139
15	B10	14-Deoxyandrographolide	31140
15	B11	Indole-3-carboxylic Acid	31166
15	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
15	C1	<i>Unused</i>	
15	C2	<i>Unused</i>	
15	C3	(±)-Epibatidine (hydrochloride)	31193
15	C4	Capsanthin	31202
15	C5	Sophoricoside	31203
15	C6	Hydroxysafflor Yellow A	31204
15	C7	4-Hydroxybenzyl Alcohol	31209
15	C8	Isovitexin	31212
15	C9	Ajmalicine	31213
15	C10	Chelidonic Acid	31324
15	C11	Nebularine	31329
15	C12	<i>Unused</i>	
15	D1	<i>Unused</i>	
15	D2	Spiculisporic Acid	31334
15	D3	Sclareolide	31371
15	D4	Aspyrone	31389
15	D5	(±)-α-Bisabolol	31397
15	D6	Anethol	31398
15	D7	Safranal	31399
15	D8	Gibberellin A ₄	31400
15	D9	Corynoxine	31403
15	D10	Ginsenoside F ₃	31409
15	D11	Acetomycin	31410
15	D12	<i>Unused</i>	
15	E1	<i>Unused</i>	
15	E2	Echinospurin	31416
15	E3	Ascochin	31450
15	E4	<i>trans</i> -β-Farnesene	31502
15	E5	Guaiazulene	31506
15	E6	Theaflavin-3-gallate	31510
15	E7	Quillaic Acid	31512
15	E8	Alyssin	31513
15	E9	Acetoxyvaleric Acid	31516
15	E10	Macranthoidin B	31533
15	E11	(+)-δ-Tocopherol	31589
15	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
15	F1	<i>Unused</i>	
15	F2	Poncirin	31706
15	F3	Delsoline	31748
15	F4	Erianin	32516
15	F5	Parvodicin C ₂	32542
15	F6	Chebulinic Acid	32554
15	F7	Dihydrofolic Acid	32750
15	F8	Sesamol	32903
15	F9	1-Hydroxyanthraquinone	32905
15	F10	2-Hydroxyanthraquinone	32906
15	F11	Se-Methylselenocysteine (hydrochloride)	32950
15	F12	<i>Unused</i>	
15	G1	<i>Unused</i>	
15	G2	(±)-Salsolidine	32953
15	G3	10-deacetyl-7-xylosyl Paclitaxel	32955
15	G4	(-)-Lupinine	33026
15	G5	<i>Unused</i>	
15	G6	<i>Unused</i>	
15	G7	<i>Unused</i>	
15	G8	28-Acetylbetulin	33052
15	G9	3-Oxobetulin	33066
15	G10	Sparstolonin B	33080
15	G11	Sophoradin	33087
15	G12	<i>Unused</i>	
15	H1	<i>Unused</i>	
15	H2	(-)-Physostigmine (salicylate)	33106
15	H3	(-)-dehydro Griseofulvin	33117
15	H4	6,2',4'-Trimethoxyflavone	33150
15	H5	(+)-Secoisolariciresinol	33155
15	H6	Saikosaponin D (hydrate)	33156
15	H7	Tetrahydroberberine	33157
15	H8	Pachymic Acid	33180
15	H9	Luteolin 7-O-Glucuronide	33188
15	H10	<i>Unused</i>	
15	H11	Rebaudioside B	33199
15	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
16	A1	<i>Unused</i>	
16	A2	20(S)-Protopanaxadiol	33201
16	A3	Asperosaponin VI	33206
16	A4	Icariside I	33209
16	A5	Clitorin	33211
16	A6	Ganoderic Acid C1	33212
16	A7	Marein	33213
16	A8	Pseudoprotodioscin	33215
16	A9	Tribuloside (hydrate)	33218
16	A10	20-Deoxyingenol	33225
16	A11	<i>Unused</i>	
16	A12	<i>Unused</i>	
16	B1	<i>Unused</i>	
16	B2	Benzoylmesaconine	33238
16	B3	Ganoderic Acid C6	33241
16	B4	<i>Unused</i>	
16	B5	Pseudomonic Acid (lithium salt)	33271
16	B6	Chrysoeriol	33395
16	B7	Aromadendrin	33402
16	B8	Vicenin-2	33403
16	B9	Nordihydrocapsaicin	33407
16	B10	Previtamin D ₃	33409
16	B11	Helicid	33414
16	B12	<i>Unused</i>	
16	C1	<i>Unused</i>	
16	C2	Kuwanon C	33421
16	C3	Asiaticoside B	33427
16	C4	<i>Unused</i>	
16	C5	(-)-Sweroside	33454
16	C6	Ginkgotoxin	33460
16	C7	Syringaldehyde	33489
16	C8	Piperitone	33490
16	C9	Dihydrocoumarin	33495
16	C10	7-Methoxyflavone	33497
16	C11	N,N'-Diphenethylurea	33580
16	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
16	D1	<i>Unused</i>	
16	D2	Tubercmycin B	33659
16	D3	FR900359	33666
16	D4	Octahydrocurcumin	33686
16	D5	Lithospermic Acid	33694
16	D6	Punicalin	33714
16	D7	Dehydroabietic Acid	33722
16	D8	Kaempferol 3-O-galactoside	33747
16	D9	Ricinine	33748
16	D10	Carvacrol	33753
16	D11	Apigenin 7-O-Glucuronide (hydrate)	33754
16	D12	<i>Unused</i>	
16	E1	<i>Unused</i>	
16	E2	Urolithin B	33757
16	E3	Myrislignan	33759
16	E4	(S)-Canadine	33805
16	E5	Naringenin Chalcone	33806
16	E6	Eupalinolide A	33808
16	E7	Perhydrofarnesyl Acetone	33825
16	E8	D-Arabinitol	33826
16	E9	Aromadendrene	33829
16	E10	Isoalantolactone	33838
16	E11	Nuciferine	33839
16	E12	<i>Unused</i>	
16	F1	<i>Unused</i>	
16	F2	Gossypetin	33840
16	F3	(+)-Oxypeucedanin hydrate	33841
16	F4	1,4-Dicaffeoylquinic Acid	33842
16	F5	Arteannuin B	33852
16	F6	Thermopsine	33853
16	F7	Scopolin	33874
16	F8	Marrubiin	33876
16	F9	Gentisein	33877
16	F10	Pyridoxamine (hydrochloride)	33899
16	F11	β -Phellandrene	33905
16	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
16	G1	<i>Unused</i>	
16	G2	(+)-Ledene	33906
16	G3	<i>Unused</i>	
16	G4	Sinapyl Alcohol	33916
16	G5	Avicularin	33918
16	G6	γ -Glu-Leu	33923
16	G7	Escin lb	33926
16	G8	Isoflavone	33964
16	G9	Tectoridin	33969
16	G10	Vitamin K ₄	33970
16	G11	Brevilin A	33993
16	G12	<i>Unused</i>	
16	H1	<i>Unused</i>	
16	H2	Flavone	34028
16	H3	Prunin	34034
16	H4	Isorhoifolin	34036
16	H5	Toosendanin	34042
16	H6	Benzoylhypaconine	34046
16	H7	Shanzhiside methyl ester	34047
16	H8	Alisol A	34050
16	H9	Escin la	34057
16	H10	Didemnin B	34063
16	H11	Xanthone	34098
16	H12	<i>Unused</i>	
17	A1	<i>Unused</i>	
17	A2	Isoscopoletin	34117
17	A3	N-Demethylvancomycin (hydrochloride)	34119
17	A4	N'-Nitrosonornicotine	34120
17	A5	<i>meta</i> -Topolin	34122
17	A6	Yunaconitine	34127
17	A7	Zingibroside R1	34130
17	A8	Visnagin	34140
17	A9	Isobergapten	34141
17	A10	Gossypin	34142
17	A11	Biotin (S)-sulfoxide	34145
17	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
17	B1	<i>Unused</i>	
17	B2	Ethyl Gallate	34193
17	B3	Loliolide	34196
17	B4	Hederacoside D	34202
17	B5	Maduramicin (ammonium salt)	34270
17	B6	Tropinone	34287
17	B7	3-Methylxanthine	34288
17	B8	Tricyclene	34290
17	B9	Gibberellin A ₁	34291
17	B10	1F-Fructofuranosylnystose	34292
17	B11	α-Cyperone	34294
17	B12	<i>Unused</i>	
17	C1	<i>Unused</i>	
17	C2	7- <i>epi</i> -10-Deacetylaxol	34296
17	C3	Notoginsenoside Fe	34297
17	C4	Ganoderol B	34324
17	C5	1-Kestose	34346
17	C6	Kaempferol 7-O-glucoside	34347
17	C7	7-Methylxanthine	34348
17	C8	5-Feruloylquinic Acid	34349
17	C9	Shikalkin	34355
17	C10	Epimedin C	34375
17	C11	Gardenoside	34381
17	C12	<i>Unused</i>	
17	D1	<i>Unused</i>	
17	D2	Quassin	34382
17	D3	Rosin	34383
17	D4	Oxyberberine	34384
17	D5	Ophiopogonin D	34385
17	D6	Levistolide A	34386
17	D7	Nepitrin	34387
17	D8	Poliumoside	34389
17	D9	Triptophenolide	34390
17	D10	Litorin (trifluoroacetate salt)	34427
17	D11	(+)-δ-Cadinene	34444
17	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
17	E1	<i>Unused</i>	
17	E2	Tuberostemonine	34457
17	E3	Bufotalin	34466
17	E4	Orcinol	34467
17	E5	8-Oxycoptisine	34472
17	E6	Emodin 8-glucoside	34474
17	E7	Tiliroside	34493
17	E8	Thymol	34496
17	E9	(±)-6-Prenylnaringenin	34499
17	E10	Gypenoside XVII	34514
17	E11	Kaempferol 3-neohesperidoside	34533
17	E12	<i>Unused</i>	
17	F1	<i>Unused</i>	
17	F2	Ganoderic Acid DM	34536
17	F3	Isopimaric Acid	34544
17	F4	Hesperetin 7-O-glucoside	34545
17	F5	<i>Unused</i>	
17	F6	Epiberberine	34569
17	F7	Byakangelicol	34571
17	F8	Caudatin	34585
17	F9	2,3,5,4'-Tetrahydroxyl-diphenylethylene 2-O-β-D-Glucoside	34594
17	F10	Apiin	34596
17	F11	Saikosaponin C	34599
17	F12	<i>Unused</i>	
17	G1	<i>Unused</i>	
17	G2	Byakangelicin	34600
17	G3	Songorine	34601
17	G4	Tetrahydrocoptisine	34603
17	G5	Isorhamnetin 3-O-neohesperidoside	34604
17	G6	Saikosaponin B ₁	34609
17	G7	Deltaline	34612
17	G8	Podocarpusflavone A	34613
17	G9	Incensole Acetate	34614
17	G10	Ardisiacrispin A	34615
17	G11	Corylifol A	34616
17	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
17	H1	<i>Unused</i>	
17	H2	Veratrosine	34617
17	H3	Isopimpinellin	34642
17	H4	Skimmianine	34643
17	H5	Kaempferol 3-glucuronide	34645
17	H6	Dictamnine	34646
17	H7	Isocorynoxine	34653
17	H8	<i>Unused</i>	
17	H9	Trifolirhizin	34655
17	H10	Siamenoside I	34656
17	H11	Engletin	34657
17	H12	<i>Unused</i>	
18	A1	<i>Unused</i>	
18	A2	Silychristin	34658
18	A3	Verbenalin	34663
18	A4	Fraxinellone	34672
18	A5	Bavachalcone	34673
18	A6	Crassicauline A	34674
18	A7	Ganoderic Acid G	34675
18	A8	Irigenin	34676
18	A9	Gitogenin	34681
18	A10	Menthone	34683
18	A11	α -Guaiene	34686
18	A12	<i>Unused</i>	
18	B1	<i>Unused</i>	
18	B2	($\pm$)-Marinopyrrole A	34723
18	B3	DL-Sulforaphane Glutathione	34731
18	B4	Meranzin	34743
18	B5	Swertisin	34744
18	B6	Disogluside	34746
18	B7	N-Methylcytisine	34752
18	B8	Neferine	34753
18	B9	Sophoraflavanone G	34755
18	B10	N-Nornuciferine	34756
18	B11	Obtusifolin	34757
18	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 47 of 59

Plate	Well	Contents	Item Number
18	C1	<i>Unused</i>	
18	C2	Cyclo(Phe-Pro)	34758
18	C3	Murrayone	34767
18	C4	Lithospermoxide	34771
18	C5	Ranaconitine	34774
18	C6	Catechol	34791
18	C7	Phellodendrine (chloride)	34802
18	C8	Decursinol	34807
18	C9	Alisol B	34808
18	C10	Blumeatin	34809
18	C11	Grosvenorine	34810
18	C12	<i>Unused</i>	
18	D1	<i>Unused</i>	
18	D2	3-Hydroxycoumarin	34815
18	D3	4-Hexylresorcinol	34824
18	D4	Chikusetsusaponin IVa	34832
18	D5	all- <i>trans</i> -13,14-Dihydroretinol	34873
18	D6	(E/Z)-Ajoene	34881
18	D7	Germacrene B	34907
18	D8	Prim-O-Glucosylcimifugin	34921
18	D9	Brevifolincarboxylic Acid	34922
18	D10	Vitexin-4''-O-glucoside	34923
18	D11	Oroxin A	34924
18	D12	<i>Unused</i>	
18	E1	<i>Unused</i>	
18	E2	Dehydrotumulosic Acid	34938
18	E3	Hirsuteine	34939
18	E4	Rubiadin-1 methyl ether	34940
18	E5	3-Butylidenephthalide	34983
18	E6	Agrimol B	34996
18	E7	Perillene	35006
18	E8	Corydine	35028
18	E9	Wilforgine	35029
18	E10	Licoflavone C	35031
18	E11	Columbamine	35032
18	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 48 of 59

Plate	Well	Contents	Item Number
18	F1	<i>Unused</i>	
18	F2	Atractylodin	35033
18	F3	Chrysin 7-Glucuronide	35034
18	F4	Suberosin	35035
18	F5	Periplogenin	35047
18	F6	Isomenthone	35063
18	F7	4-Hydroxyquinoline	35064
18	F8	N-desmethyl Galantamine	35074
18	F9	O-desmethyl Galantamine	35080
18	F10	Pyrrolnitrin	35099
18	F11	Lotusine	35136
18	F12	<i>Unused</i>	
18	G1	<i>Unused</i>	
18	G2	1-O-Acetylbritannilactone	35137
18	G3	(2S)-6-Prenylnaringenin	35138
18	G4	Scoulerine	35140
18	G5	Quercetin 3-O-β-D-xylopyranoside	35141
18	G6	Senkyunolide H	35142
18	G7	Rosmanol	35143
18	G8	Akuammine	35149
18	G9	Methyl 4-hydroxycinnamate	35203
18	G10	Rosarin	35214
18	G11	Hypaphorine	35228
18	G12	<i>Unused</i>	
18	H1	<i>Unused</i>	
18	H2	Genkwanin	35229
18	H3	5-O-Demethylnobiletin	35230
18	H4	N-methyl Tyramine (hydrochloride)	35233
18	H5	Sennidin A	35237
18	H6	Rebaudioside M	35239
18	H7	Continentalic Acid	35240
18	H8	1-Methylxanthine	35245
18	H9	Galantamine (hydrobromide)	35247
18	H10	Thujopsene	35270
18	H11	Robinetin	35272
18	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 49 of 59

Plate	Well	Contents	Item Number
19	A1	<i>Unused</i>	
19	A2	7-Demethylsuberosin	35282
19	A3	Scopoline	35301
19	A4	Ginkgolide K	35304
19	A5	Larixyl Acetate	35309
19	A6	Lupenone	35335
19	A7	N ⁶ ,N ⁶ -Dimethyladenosine	35348
19	A8	Dihydrojasmonic Acid	35362
19	A9	Paederosidic Acid	35363
19	A10	α-Arbutin	35386
19	A11	Xylotetraose	35387
19	A12	<i>Unused</i>	
19	B1	<i>Unused</i>	
19	B2	7,4'-Di-O-methylapigenin	35388
19	B3	1-Methylpseudouridine	35401
19	B4	6,7,4'-Trihydroxyisoflavone	35405
19	B5	Dihydronarwedine	35409
19	B6	Galanthamine N-oxide	35413
19	B7	Indican	35417
19	B8	Galgravin	35431
19	B9	Primin	35441
19	B10	7,4'-Dihydroxyflavone	35491
19	B11	Aristolochic Acid D	35492
19	B12	<i>Unused</i>	
19	C1	<i>Unused</i>	
19	C2	20(R)-Ginsenoside Rh ₁	35493
19	C3	Kumatakenin	35514
19	C4	4-Chloroindole-3-acetic Acid	35534
19	C5	Oleuropein aglycone	35547
19	C6	cis-Zeatin	35549
19	C7	1-Octen-3-ol	35550
19	C8	Hygromycin A	35565
19	C9	D-Tetrahydropalmatine	35577
19	C10	Rhoifolin	35583
19	C11	Berbamine (hydrochloride)	35593
19	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
19	D1	<i>Unused</i>	
19	D2	1-Isothiocyanato-6-(methylsulphenyl)-hexane	35595
19	D3	Isorhamnetin-3-O- β -D-Glucoside	35600
19	D4	Epigoitrin	35620
19	D5	Genipin 1-gentiobioside	35621
19	D6	2-keto D-Gluconic Acid (calcium salt)	35622
19	D7	Tirucallol	35623
19	D8	Dipsacoside B	35624
19	D9	Saikosaponin F	35626
19	D10	Alisol F	35627
19	D11	Camelliaside A	35628
19	D12	<i>Unused</i>	
19	E1	<i>Unused</i>	
19	E2	Huperzine B	35629
19	E3	Gartanin	35631
19	E4	Salvigenin	35632
19	E5	4-Acetylphenol	35636
19	E6	Senkyunolide I	35649
19	E7	Coclaurine	35716
19	E8	Ginkgolic Acid C17:1	35737
19	E9	($\pm$)-Tuberonic Acid	35750
19	E10	<i>Unused</i>	
19	E11	Techtochrysin	35791
19	E12	<i>Unused</i>	
19	F1	<i>Unused</i>	
19	F2	Saponarin	35805
19	F3	Ononetin	35813
19	F4	4-Methoxyphenyl- β -D-Glucopyranoside	35818
19	F5	4-oxo Retinoic Acid	35819
19	F6	Bayogenin	35822
19	F7	Alisol C Monoacetate	35823
19	F8	Isomeranzin	35827
19	F9	N-acetyl-2-Aminophenol	35829
19	F10	Tubercidin	35834
19	F11	Vanillin 4-O- β -D-Glucoside	35908
19	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 51 of 59

Plate	Well	Contents	Item Number
19	G1	<i>Unused</i>	
19	G2	Eriodictyol 7-O-glucoside	36003
19	G3	Yamogenin	36008
19	G4	Isosinensetin	36009
19	G5	β -Glucogallin	36010
19	G6	Cycleanine	36020
19	G7	Isolariciresinol	36032
19	G8	Remerin	36035
19	G9	α -Cembrenediol	36042
19	G10	7-Hydroxyflavone	36071
19	G11	Chamazulene	36072
19	G12	<i>Unused</i>	
19	H1	<i>Unused</i>	
19	H2	Flavokawain A	36083
19	H3	Frangulin A	36087
19	H4	Harmalol (hydrochloride hydrate)	36090
19	H5	Quercetin 3-O- β -D-Glucose-7-O- β -D-Gentiobioside	36130
19	H6	<i>Unused</i>	
19	H7	<i>Unused</i>	
19	H8	3'-Hydroxygenkwanin	36150
19	H9	3-Deoxyaconitine	36151
19	H10	3-hydroxy-9,10-Dimethoxypterocarpan	36152
19	H11	3-Acetylaconitine	36155
19	H12	<i>Unused</i>	
20	A1	<i>Unused</i>	
20	A2	2"-O- β -L-Galactopyranosylorientin	36156
20	A3	Aristolone	36157
20	A4	3,29-Dibenzoyl Rarounitriol	36158
20	A5	Euscaphic Acid	36162
20	A6	1,3-Dimethyluric Acid	36167
20	A7	Bullatine B	36170
20	A8	10-Shogaol	36171
20	A9	Cucurbitacin D	36172
20	A10	Atractylenolide II	36173
20	A11	Timosaponin AI	36174
20	A12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 52 of 59

Plate	Well	Contents	Item Number
20	B1	<i>Unused</i>	
20	B2	12-O-Methylcarnosic Acid	36175
20	B3	Crebanine	36176
20	B4	<i>Unused</i>	
20	B5	γ -Glu-Phe (trifluoroacetate salt)	36181
20	B6	1-O-Caffeoylquinic Acid	36189
20	B7	4',7-Di-O-methylnaringenin	36190
20	B8	Androsin	36191
20	B9	9,10-Dimethoxypterocarpan-3-O- β -D-glucoside	36192
20	B10	Arnicolide D	36193
20	B11	Clinodiside A	36200
20	B12	<i>Unused</i>	
20	C1	<i>Unused</i>	
20	C2	Koumine	36210
20	C3	D-Mannosamine (hydrochloride)	36231
20	C4	3-(3-Hydroxyphenyl)propanoic Acid	36268
20	C5	Rhamnocitrin	36272
20	C6	Fustin	36273
20	C7	Pimpinellin	36274
20	C8	Cornuside	36284
20	C9	Parishin B	36296
20	C10	Hinokiflavone	36338
20	C11	Pulmatin	36343
20	C12	<i>Unused</i>	
20	D1	<i>Unused</i>	
20	D2	Methylprotodioscin	36344
20	D3	Rhodojaponin III	36345
20	D4	(+)- γ -Tocopherol	36362
20	D5	3-(4-Hydroxyphenyl)lactic Acid	36367
20	D6	DL-Pyroglutamic Acid	36375
20	D7	Hexahydrohippuric Acid	36386
20	D8	Procyanidin B4	36398
20	D9	(+)-Sparteine	36400
20	D10	2-Methoxy-4-vinylphenol	36427
20	D11	4-Methylcatechol	36429
20	D12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 53 of 59

Plate	Well	Contents	Item Number
20	E1	<i>Unused</i>	
20	E2	Deoxypodophyllotoxin	36434
20	E3	O-Desmethylangolensin	36435
20	E4	5,6-Dimethylbenzimidazole	36437
20	E5	Myricetin 3-galactoside	36467
20	E6	3,5,6,7,8,3',4'-Heptamethoxyflavone	36475
20	E7	4'-Demethylpodophyllotoxin	36481
20	E8	1-Cinnamoylpyrrolidine	36483
20	E9	3-Isomangostin	36484
20	E10	Dimethylfraxetin	36485
20	E11	2'-Acetylacteoside	36511
20	E12	<i>Unused</i>	
20	F1	<i>Unused</i>	
20	F2	Cephalotaxine	36521
20	F3	Valtrate	36522
20	F4	Neoglycyrol	36525
20	F5	3- <i>epi</i> Ursolic Acid	36528
20	F6	4-Allylcatechol	36542
20	F7	12-Epinapelline	36550
20	F8	Sulfuretin	36569
20	F9	1,1,1,1-Kestohexose	36591
20	F10	Miltirone	36663
20	F11	(-)-Syringaresinol di-O-glucoside	36672
20	F12	<i>Unused</i>	
20	G1	<i>Unused</i>	
20	G2	γ -Glu-Ala (trifluoroacetate salt)	36688
20	G3	Aposcopolamine	36689
20	G4	5-(3',4'-Dihydroxyphenyl)- γ -Valerolactone	36690
20	G5	<i>Unused</i>	
20	G6	Brazilein	36736
20	G7	3'-hydroxy Puerarin	36771
20	G8	Butyl 4-hydroxybenzoate	36797
20	G9	Asaraldehyde	36798
20	G10	N-formyl-2-Aminophenol	36802
20	G11	Gingerenone A	36841
20	G12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 54 of 59

Plate	Well	Contents	Item Number
20	H1	<i>Unused</i>	
20	H2	Febrifugine (hydrochloride)	37044
20	H3	<i>Unused</i>	
20	H4	Ziyuglycoside I	37320
20	H5	Ethylparaben	37357
20	H6	Hispidol	37373
20	H7	<i>Unused</i>	
20	H8	Propylparaben	37417
20	H9	Ermanin	37535
20	H10	Lentztrehalose A	37732
20	H11	Lentztrehalose B	37733
20	H12	<i>Unused</i>	
21	A1	<i>Unused</i>	
21	A2	Rhyncophylline	37946
21	A3	Mycophenolic Acid methyl ester	38397
21	A4	5-(3'-Hydroxyphenyl)- γ -Valerolactone	38519
21	A5	Bergaptol	38742
21	A6	Myriocin	63150
21	A7	<i>Unused</i>	
21	A8	Parthenolide	70080
21	A9	Guaiacol	70430
21	A10	Filipin III	70440
21	A11	Bestatin (hydrochloride)	70520
21	A12	<i>Unused</i>	
21	B1	<i>Unused</i>	
21	B2	3,4-Dihydroxyphenyl ethanol	70604
21	B3	Baicalein	70610
21	B4	<i>trans</i> -Resveratrol	70675
21	B5	<i>rac</i> -Astaxanthin	70685
21	B6	Rosmarinic Acid	70900
21	B7	Chlorogenic Acid	70930
21	B8	(-)-Epigallocatechin Gallate	70935
21	B9	(+)-Catechin (hydrate)	70940
21	B10	<i>Unused</i>	
21	B11	Celastrol	70950
21	B12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Plate	Well	Contents	Item Number
21	C1	<i>Unused</i>	
21	C2	Lactacystin	70980
21	C3	(Z)-Guggulsterone	71800
21	C4	<i>Unused</i>	
21	C5	Curcumin	81025
21	C6	(±)-Jasmonic Acid	88300
21	C7	γ-CEHC	89630
21	C8	Trichostatin A	89730
21	C9	Carnosol	89800
21	C10	Carnosic Acid	89820
21	C11	Capsaicin	92350
21	C12	<i>Unused</i>	
21	D1	<i>Unused</i>	
21	D2	Dihydrocapsaicin	92355
21	D3	8-methyl Nonanoic Acid	9000310
21	D4	Pepstatin A	9000469
21	D5	25-hydroxy Vitamin D ₃	9000683
21	D6	4-chloro-α-Cyanocinnamic Acid	9000981
21	D7	N-3-hydroxydecanoyl-L-Homoserine lactone	9001147
21	D8	Piperlonguminine	9001155
21	D9	Resibufogenin	9001348
21	D10	Nanaomycin A	9001479
21	D11	β-Tocotrienol	9001533
21	D12	<i>Unused</i>	
21	E1	<i>Unused</i>	
21	E2	Salvianolic Acid B	9001577
21	E3	Propionyl-L-carnitine (chloride)	9001873
21	E4	(-)-α-Bisabolol	9002495
21	E5	<i>Unused</i>	
21	E6	Geranylgeranoic Acid	9003474
21	E7	2-Nonylquinolin-4(1H)-one	9003627
21	E8	L-3-n-Butylphthalide	9003759
21	E9	3-(3-Hydroxyphenyl)propionic Acid sulfate	9004103
21	E10	p-Carboxyphenyl Sulfate	9004155
21	E11	Luteolin	10004161
21	E12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 56 of 59

Plate	Well	Contents	Item Number
21	F1	<i>Unused</i>	
21	F2	3,4',5-Trimethoxybenzophenone	10004185
21	F3	cis-Resveratrol	10004235
21	F4	Ionomycin	10004974
21	F5	Daidzein	10005166
21	F6	Genistein	10005167
21	F7	Quercetin (hydrate)	10005169
21	F8	Matairesinol	10005174
21	F9	Cerulenin	10005647
21	F10	Hesperetin	10006084
21	F11	Leelamine	10006148
21	F12	<i>Unused</i>	
21	G1	<i>Unused</i>	
21	G2	Silybin	10006211
21	G3	Berberine (chloride)	10006427
21	G4	E-64	10007963
21	G5	FK-506	10007965
21	G6	α -Tocotrienol	10008377
21	G7	γ -Tocotrienol	10008494
21	G8	δ -Tocotrienol	10008513
21	G9	Leelamine (hydrochloride)	10008614
21	G10	<i>Unused</i>	
21	G11	(+)-2,5- <i>epi</i> Goniethalesdiol	10008887
21	G12	<i>Unused</i>	
21	H1	<i>Unused</i>	
21	H2	Piceatannol	10009366
21	H3	Kifunensine	10009437
21	H4	Bergamottin	10009439
21	H5	<i>Unused</i>	
21	H6	Pyocyanin	10009594
21	H7	Maslinic Acid	10009645
21	H8	Tangeretin	10009911
21	H9	Zeaxanthin	10009992
21	H10	Inotilone	10010089
21	H11	(R)-Equol	10010172
21	H12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 57 of 59

Plate	Well	Contents	Item Number
22	A1	<i>Unused</i>	
22	A2	(S)-Equol	10010173
22	A3	Lavendustin A	10010268
22	A4	Apigenin	10010275
22	A5	Harmine	10010324
22	A6	Lovastatin	10010338
22	A7	Lovastatin Hydroxy Acid (sodium salt)	10010339
22	A8	Vialinin A	10010519
22	A9	Genipin	10010622
22	A10	(-)-Deguelin	10010706
22	A11	Lutein	10010811
22	A12	<i>Unused</i>	
22	B1	<i>Unused</i>	
22	B2	Steviol	10011344
22	B3	1-Deoxynojirimycin (hydrochloride)	10011718
22	B4	Myricetin	10012600
22	B5	Hispidin	10012605
22	B6	Cyclosporin H	17182
22	B7	<i>Unused</i>	
22	B8	<i>Unused</i>	
22	B9	<i>Unused</i>	
22	B10	<i>Unused</i>	
22	B11	<i>Unused</i>	
22	B12	<i>Unused</i>	
22	C1	<i>Unused</i>	
22	C2	<i>Unused</i>	
22	C3	<i>Unused</i>	
22	C4	<i>Unused</i>	
22	C5	<i>Unused</i>	
22	C6	<i>Unused</i>	
22	C7	<i>Unused</i>	
22	C8	<i>Unused</i>	
22	C9	<i>Unused</i>	
22	C10	<i>Unused</i>	
22	C11	<i>Unused</i>	
22	C12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Page 58 of 59

Plate	Well	Contents	Item Number
22	D1	<i>Unused</i>	
22	D2	Oligomycin Complex	11341
22	D3	Concanavalin A	14951
22	D4	Moenomycin Complex	15506
22	D5	Teicoplanin Complex	15511
22	D6	Bromelain	17337
22	D7	Carotenoid Mixture	18264
22	D8	Bacitracin (zinc)	21212
22	D9	Tyrothricin	22177
22	D10	Tylosin	22599
22	D11	Crocin/Gardenia Extract	26337
22	D12	<i>Unused</i>	
22	E1	<i>Unused</i>	
22	E2	Ramoplanin Complex	28189
22	E3	Hesperidin methylchalcone/ <i>C. aurantium</i> L. Extract (hydrate)	30884
22	E4	Antimycin A Complex	34799
22	E5	Enniatin Complex	9002040
22	E6	<i>Unused</i>	
22	E7	<i>Unused</i>	
22	E8	<i>Unused</i>	
22	E9	<i>Unused</i>	
22	E10	<i>Unused</i>	
22	E11	<i>Unused</i>	
22	E12	<i>Unused</i>	
22	F1	<i>Unused</i>	
22	F2	<i>Unused</i>	
22	F3	<i>Unused</i>	
22	F4	<i>Unused</i>	
22	F5	<i>Unused</i>	
22	F6	<i>Unused</i>	
22	F7	<i>Unused</i>	
22	F8	<i>Unused</i>	
22	F9	<i>Unused</i>	
22	F10	<i>Unused</i>	
22	F11	<i>Unused</i>	
22	F12	<i>Unused</i>	

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

[illegible]

WWW.CAYMANCHEM.COM