

PRODUCT INFORMATION



Bio-Active Lipid I Screening Library (96-Well)

Item No. 10506 • Batch No. 0608063

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

Page 1 of 35

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	Prostaglandin A ₁	10010
1	A3	1-Arachidonoyl Lysophosphatidic Acid (sodium salt)	10019
1	A4	POV-PC	10031
1	A5	N-(α -Linolenoyl) Tyrosine	10032
1	A6	10-Nitrolinoleate	10037
1	A7	PGPC	10044
1	A8	15-deoxy- $\Delta^{12,14}$ -Prostaglandin A ₁	10065
1	A9	15- <i>epi</i> Prostaglandin A ₁	10070
1	A10	16,16-dimethyl Prostaglandin A ₁	10080
1	A11	N-Oleoyl Dopamine	10115
1	A12	Unused	
1	B1	Unused	
1	B2	Prostaglandin F _{2α} -1-glycerol ester	10139
1	B3	Prostaglandin E ₂ -1-glycerol ester	10140
1	B4	N-Arachidonoyl-3-hydroxy- γ -Aminobutyric Acid	10158
1	B5	Arachidonoyl <i>p</i> -Nitroaniline	10168
1	B6	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PC	10172
1	B7	Arachidonoyl Ethanolamide Phosphate	10180
1	B8	Prostaglandin D ₂ serinol amide	10192
1	B9	Prostaglandin E ₂ serinol amide	10193
1	B10	Prostaglandin F _{2α} serinol amide	10194
1	B11	Prostaglandin A ₂	10210
1	B12	Unused	
1	C1	Unused	
1	C2	13,14-dihydro-15-keto Prostaglandin A ₂	10260
1	C3	15-deoxy- $\Delta^{12,14}$ -Prostaglandin A ₂	10265
1	C4	16,16-dimethyl Prostaglandin A ₂	10280
1	C5	16-phenoxy tetranor Prostaglandin A ₂	10285
1	C6	17-phenyl trinor Prostaglandin A ₂	10288
1	C7	17-phenyl trinor-13,14-dihydro Prostaglandin A ₂	10290
1	C8	7-hydroxycoumarinyl- γ -Linolenate	10556
1	C9	all- <i>cis</i> -14,17,20,23,26,29-Dotriacontahexaenol	10562
1	C10	Heneicosapentaenoic Acid	10670
1	C11	($\pm$)-Jasmonic Acid-Isoleucine	10740
1	C12	Unused	

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, 06/25/2025

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 35

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Eicosapentaenoyl Ethanolamide	10964
1	D3	Palmitoleoyl Ethanolamide	10965
1	D4	10-HDA	10976
1	D5	Nocloprost	10988
1	D6	18-methyl Eicosanoic Acid	11009
1	D7	Palmitoyl-DL-carnitine (chloride)	11095
1	D8	25-hydroxy Cholesterol	11097
1	D9	Prostaglandin B ₁	11110
1	D10	Δ^2 - <i>trans</i> -Hexadecenoic Acid	11132
1	D11	Δ^2 - <i>cis</i> -Hexadecenoic Acid	11133
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Termitomycamide B	11190
1	E3	Termitomycamide E	11191
1	E4	Prostaglandin B ₂	11210
1	E5	Sulfosuccinimidyl Oleate (sodium salt)	11211
1	E6	Tafluprost ethyl ester	11612
1	E7	11-keto- β -Boswellic Acid	11668
1	E8	3-acetyl-11-keto- β -Boswellic Acid	11672
1	E9	β -Boswellic Acid	11693
1	E10	24 α -ethyl Cholesterol (technical grade)	11756
1	E11	L- α -Hydroxyglutaric Acid (sodium salt)	11876
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	2-O-Hexadecyl Glycerol	11924
1	F3	Treprostinil (diethanolamine salt)	11927
1	F4	<i>cis</i> -2-Decenoic Acid	11966
1	F5	Octyl- α -ketoglutarate	11970
1	F6	Prostaglandin D ₁	12000
1	F7	Prostaglandin D ₁ Alcohol	12002
1	F8	Prostaglandin D ₂	12010
1	F9	Prostaglandin D ₂ Ethanolamide	12012
1	F10	11-deoxy-11-methylene Prostaglandin D ₂	12410
1	F11	13,14-dihydro-15-keto Prostaglandin D ₂	12610
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



Page 3 of 35

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	15-deoxy- $\Delta^{12,14}$ -Prostaglandin D ₂	12700
1	G3	15(S)-15-methyl Prostaglandin D ₂	12730
1	G4	16,16-dimethyl Prostaglandin D ₂	12750
1	G5	17-phenyl trinor Prostaglandin D ₂	12810
1	G6	Prostaglandin E ₁ Ethanolamide	13012
1	G7	2-Hydroxyestradiol	13019
1	G8	Prostaglandin E ₁ Alcohol	13020
1	G9	2-Methoxyestradiol	13021
1	G10	1a,1b-dihomo Prostaglandin E ₁	13050
1	G11	N-Oleoyl-L-Serine	13058
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Anacardic Acid	13144
1	H3	9(Z)-Tricosene	13236
1	H4	Farnesyl Alcohol	13268
1	H5	Myristic Acid	13351
1	H6	Lignoceric Acid	13353
1	H7	Hexacosanoic Acid	13354
1	H8	8-iso Prostaglandin E ₁	13360
1	H9	11 β -Prostaglandin E ₁	13450
1	H10	11-deoxy Prostaglandin E ₁	13510
1	H11	1-Deoxysphinganine (m18:0)	13511
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	17-phenyl trinor Prostaglandin E ₂ ethyl amide	13532
2	A3	13,14-dihydro-16,16-difluoro Prostaglandin F _{2α}	13609
2	A4	13,14-dihydro Prostaglandin E ₁	13610
2	A5	13,14-dihydro-16,16-difluoro Prostaglandin D ₂	13612
2	A6	13,14-dihydro-15(R)-Prostaglandin E ₁	13630
2	A7	13,14-dihydro-16,16-difluoro Prostaglandin J ₂	13635
2	A8	Oleoyl Serinol	13637
2	A9	13,14-dihydro-15-keto Prostaglandin E ₁	13650
2	A10	13,14-epoxy Fluprostenol isopropyl ester	13679
2	A11	15-keto Prostaglandin E ₁	13680
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 35

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	15(S)-15-methyl Prostaglandin E ₁	13730
2	B3	16,16-dimethyl Prostaglandin E ₁	13750
2	B4	Misoprostol	13820
2	B5	Nervonic Acid	13940
2	B6	Prostaglandin E ₂	14010
2	B7	Prostaglandin E ₂ Ethanolamide	14012
2	B8	(25S)- Δ^7 -Dafachronic Acid	14101
2	B9	Trilostane	14164
2	B10	8-iso Prostaglandin E ₂	14350
2	B11	9-deoxy-9-methylene Prostaglandin E ₂	14410
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	11-deoxy Prostaglandin E ₂	14520
2	C3	11-deoxy-16,16-dimethyl Prostaglandin E ₂	14570
2	C4	7-dehydro Cholesterol	14612
2	C5	13,14-dihydro-15-keto Prostaglandin E ₂	14650
2	C6	1',4"-Sophorolactone 6',6"-diacetate	14718
2	C7	15-keto Prostaglandin E ₂	14720
2	C8	Docosapentaenoyl Ethanolamide	14724
2	C9	16,16-dimethyl Prostaglandin E ₂	14750
2	C10	16-phenoxy tetranor Prostaglandin E ₂	14760
2	C11	25(R)-27-hydroxy Cholesterol	14790
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	25(S)-27-hydroxy Cholesterol	14791
2	D3	3-oxo Stearic Acid	14807
2	D4	17-phenyl trinor Prostaglandin E ₂	14810
2	D5	Estriol 3- β -D-Glucuronide (sodium salt)	14827
2	D6	19(R)-hydroxy Prostaglandin E ₂	14910
2	D7	Finasteride	14938
2	D8	20-ethyl Prostaglandin E ₂	14940
2	D9	Exemestane	15008
2	D10	Prostaglandin F _{1α}	15010
2	D11	Prostaglandin F _{1α} Alcohol	15020
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



Page 5 of 35

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	Octanoyl-DL-carnitine (chloride)	15048
2	E3	Cholesterol Sulfate (sodium salt)	15106
2	E4	1-Stearoyl-2-hydroxy-sn-glycero-3-PC	15107
2	E5	Ursodeoxycholic Acid (sodium salt)	15121
2	E6	3-methoxy Prostaglandin F _{1α}	15170
2	E7	Pentacosanoic Acid	15197
2	E8	19-hydroxy Cholesterol	15209
2	E9	6-keto Prostaglandin F _{1α}	15210
2	E10	6,15-diketo-13,14-dihydro Prostaglandin F _{1α}	15270
2	E11	Aldosterone	15273
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	<i>trans</i> -Vaccenic Acid	15301
2	F3	Ginsenoside Rg ₁	15315
2	F4	Ginsenoside Rb ₁	15319
2	F5	Ginsenoside Rd	15329
2	F6	Ginsenoside Re	15330
2	F7	Ginsenoside Rg ₂	15331
2	F8	20(2)-Ginsenoside Rg ₃	15332
2	F9	PAF C-16 Carboxylic Acid	15392
2	F10	Prostaglandin F _{1β}	15410
2	F11	11β-Prostaglandin F _{1β}	15450
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	11-deoxy Prostaglandin F _{1β}	15510
2	G3	11-deoxy Prostaglandin F _{1α}	15512
2	G4	Ganglioside G _{T1b} Mixture (sodium salt)	15588
2	G5	13,14-dihydro Prostaglandin F _{1α}	15610
2	G6	13,14-dihydro-15-keto Prostaglandin F _{1α}	15670
2	G7	15-keto Prostaglandin F _{1α}	15710
2	G8	Bufalin	15725
2	G9	Dehydroepiandrosterone	15728
2	G10	17,20-dimethyl Prostaglandin F _{1α}	15790
2	G11	Androsterone	15872
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 35

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	Dehydroepiandrosterone Sulfate (sodium salt)	15873
2	H3	Progesterone	15876
2	H4	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{1α}	15895
2	H5	Taurodeoxycholic Acid (sodium salt hydrate)	15935
2	H6	Lauroyl-DL-Carnitine (chloride)	16006
2	H7	Prostaglandin F _{2α}	16010
2	H8	Prostaglandin F _{2α} Alcohol	16012
2	H9	Prostaglandin F _{2α} Ethanolamide	16013
2	H10	Prostaglandin F _{2α} ethyl amide	16016
2	H11	Prostaglandin F _{2α} diethyl amide	16023
2	H12	<i>Unused</i>	
3	A1	<i>Unused</i>	
3	A2	Jacaric Acid	16036
3	A3	Corticosterone	16063
3	A4	20-hydroxy Ecdysone	16145
3	A5	17α,20β-Dihydroxy-4-pregnen-3-one	16146
3	A6	Estradiol 3-(β-D-Glucuronide) (sodium salt)	16155
3	A7	Estradiol 17-(β-D-Glucuronide) (sodium salt hydrate)	16156
3	A8	Taurocholic Acid (sodium salt)	16215
3	A9	7-keto Cholesterol	16339
3	A10	8-iso Prostaglandin F _{2α}	16350
3	A11	8-iso Prostaglandin F _{2β}	16370
3	A12	<i>Unused</i>	
3	B1	<i>Unused</i>	
3	B2	α-Hydroxyglutaric Acid (sodium salt)	16374
3	B3	Ponasterone A	16386
3	B4	Prostaglandin F _{2β}	16410
3	B5	16,16-dimethyl Prostaglandin F _{2β}	16470
3	B6	11-deoxy Prostaglandin F _{2α}	16500
3	B7	11-deoxy Prostaglandin F _{2β}	16510
3	B8	6-Azidohexanoic Acid	16518
3	B9	11β-Prostaglandin F _{2α}	16520
3	B10	2-Oleoyl Glycerol	16537
3	B11	Norgestimate	16547
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 35

Plate	Well	Contents	Item Number
3	C1	<i>Unused</i>	
3	C2	<i>cis,cis</i> -Octadeca-9,12-dienol	16556
3	C3	13,14-dihydro Prostaglandin F _{2α}	16660
3	C4	13,14-dihydro-15-keto Prostaglandin F _{2α}	16670
3	C5	Unoprostone	16680
3	C6	15-keto Prostaglandin F _{2α}	16720
3	C7	15(S)-15-methyl Prostaglandin F _{2α}	16743
3	C8	15(S)-15-methyl Prostaglandin F _{2α} methyl ester	16744
3	C9	16,16-dimethyl Prostaglandin F _{2α}	16750
3	C10	16-phenoxy tetranor Prostaglandin F _{2α}	16760
3	C11	(+)-Cloprostenol	16765
3	C12	<i>Unused</i>	
3	D1	<i>Unused</i>	
3	D2	Fluprostenol isopropyl ester	16769
3	D3	9-keto Fluprostenol	16781
3	D4	11-keto Fluprostenol	16783
3	D5	17-phenyl trinor Prostaglandin F _{2α}	16810
3	D6	Latanoprost (free acid)	16811
3	D7	Latanoprost	16812
3	D8	17-phenyl trinor Prostaglandin F _{2α} ethyl amide	16820
3	D9	17-phenyl trinor Prostaglandin F _{2α} amide	16821
3	D10	Latanoprost ethyl amide	16822
3	D11	17-phenyl trinor Prostaglandin F _{2α} diethyl amide	16823
3	D12	<i>Unused</i>	
3	E1	<i>Unused</i>	
3	E2	17-phenyl trinor Prostaglandin F _{2α} isopropyl ester	16824
3	E3	CAY10580	16835
3	E4	β-Carotene	16837
3	E5	13,14-dihydro-15-keto-tetranor Prostaglandin F _{1β}	16853
3	E6	(R)-3-hydroxy Myristic Acid	16871
3	E7	17-trifluoromethylphenyl trinor Prostaglandin F _{2α}	16890
3	E8	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{2α}	16895
3	E9	20-ethyl Prostaglandin F _{2α}	16940
3	E10	Glycochenodeoxycholic Acid (sodium salt hydrate)	16942
3	E11	DC-Chol (hydrochloride)	16943
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 35

Plate	Well	Contents	Item Number
3	F1	<i>Unused</i>	
3	F2	20-hydroxy Prostaglandin F _{2α}	16950
3	F3	9-PAHSA	17037
3	F4	9-PAHPA	17038
3	F5	9-OAHSA	17039
3	F6	9-POHSA	17040
3	F7	5-PAHSA	17043
3	F8	13-PAHSA	17044
3	F9	12-POHSA	17105
3	F10	12-SAHSA	17106
3	F11	12-PAHSA	17107
3	F12	<i>Unused</i>	
3	G1	<i>Unused</i>	
3	G2	12-OAHSA	17108
3	G3	9-SAHSA	17109
3	G4	13-SAHSA	17110
3	G5	13-POHSA	17111
3	G6	13-OAHSA	17112
3	G7	5-SAHSA	17113
3	G8	5-OAHSA	17115
3	G9	1-Linoleoyl-2-Hydroxy- <i>sn</i> -glycero-3-PC	17153
3	G10	Taurolithocholic Acid (sodium salt)	17275
3	G11	24-methyl Cholesterol	17344
3	G12	<i>Unused</i>	
3	H1	<i>Unused</i>	
3	H2	Pentadecanoic Acid	17399
3	H3	Isobutyryl-L-carnitine	17440
3	H4	Pyrrolidine Linoleamide	17539
3	H5	(S)-α-Methylbenzyl Ricinoleamide	17546
3	H6	Eicosapentaenoyl 1-propanol-2-amide	17547
3	H7	Pyrrolidine Ricinoleamide	17548
3	H8	(E)-2-Hexadecenal	17566
3	H9	2-Palmitoyl Glycerol	17882
3	H10	<i>trans</i> -2-Decenoic Acid	17888
3	H11	9(S)-PAHSA	18023
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	Triamcinolone acetonide	18026
4	A3	Stigmasterol	18079
4	A4	6 α -Prostaglandin I ₁	18110
4	A5	6 β -Prostaglandin I ₁	18120
4	A6	5-cis Carbaprostacyclin	18211
4	A7	13,14-dehydro-15-cyclohexyl Carbaprostacyclin	18212
4	A8	Iloprost	18215
4	A9	Ciprostene (calcium salt)	18216
4	A10	Prostaglandin I ₂ (sodium salt)	18220
4	A11	Hydrocortisone 21-hemisuccinate (sodium salt)	18226
4	A12	<i>Unused</i>	
4	B1	<i>Unused</i>	
4	B2	Ciclesonide	18227
4	B3	Beraprost (sodium salt)	18230
4	B4	Estetrol (hydrate)	18232
4	B5	5 α -Androst-16-en-3-one	18233
4	B6	Prostaglandin I ₃ (sodium salt)	18300
4	B7	($\pm$)5(6)-DiHETE lactone	18343
4	B8	($\pm$)4(5)-DiHDPA lactone	18344
4	B9	Commendamide	18421
4	B10	Anacardic Acid monoene	18422
4	B11	3 β ,5 α -THDOC	18431
4	B12	<i>Unused</i>	
4	C1	<i>Unused</i>	
4	C2	Taurolithocholic Acid 3-sulfate (sodium salt)	18468
4	C3	Prostaglandin J ₂	18500
4	C4	Δ^{12} -Prostaglandin J ₂	18550
4	C5	15-deoxy- $\Delta^{12,14}$ -Prostaglandin J ₂	18570
4	C6	CAY10410	18590
4	C7	1a,1b-dihomo Prostaglandin E ₂	18665
4	C8	PAF C-18:1	18779
4	C9	Prostaglandin K ₁	18800
4	C10	Prostaglandin K ₂	18900
4	C11	Pinane Thromboxane A ₂	19020
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	Thromboxane B ₂	19030
4	D3	12-doxy Stearic Acid	19259
4	D4	Icosabutate	19386
4	D5	11-dehydro Thromboxane B ₂	19500
4	D6	4β-hydroxy Cholesterol	19518
4	D7	6α-methyl-11β-hydroxy Progesterone	19533
4	D8	C18 Ceramide (d18:1/18:0)	19556
4	D9	(5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic Acid	19557
4	D10	Ganglioside G _{M1} (porcine brain) (sodium salt)	19579
4	D11	MMP-2 Inhibitor I	19644
4	D12	<i>Unused</i>	
4	E1	<i>Unused</i>	
4	E2	Dehydroergosterol	19709
4	E3	Undecanoic Acid	19721
4	E4	Heptadecanoic Acid	19722
4	E5	Nonadecanoic Acid	19723
4	E6	14-Pentadecenoic Acid	19724
4	E7	Tridecanoic Acid	19725
4	E8	cis-10-Heptadecenoic Acid	19748
4	E9	cis-10-Nonadecenoic Acid	19749
4	E10	Cinobufagin	19844
4	E11	Ergosterol	19850
4	E12	<i>Unused</i>	
4	F1	<i>Unused</i>	
4	F2	Pregnenolone	19864
4	F3	3-hydroxy Palmitic Acid	19934
4	F4	N-Palmitoyl-L-Aspartate	19952
4	F5	10-PAHSA	19973
4	F6	10-POHSA	19974
4	F7	10-OAHSA	19975
4	F8	10-SAHSA	19976
4	F9	cis-Vaccenic Acid	20023
4	F10	cis-Petroselinic Acid	20024
4	F11	trans-Petroselinic Acid	20026
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



Page 11 of 35

Plate	Well	Contents	Item Number
4	G1	<i>Unused</i>	
4	G2	Linoleoyl Phenylalanine	20063
4	G3	N-Oleoyl Leucine	20064
4	G4	N-Oleoyl Valine	20065
4	G5	7 α -hydroxy Cholesterol	20098
4	G6	7 β -hydroxy Cholesterol	20099
4	G7	7 β ,27-dihydroxy Cholesterol	20102
4	G8	20(S)-hydroxy Cholesterol	20103
4	G9	NSC 12	20117
4	G10	C2 dihydro Ceramide (d18:0/2:0)	20168
4	G11	Phytosphingosine (yeast)	20217
4	G12	<i>Unused</i>	
4	H1	<i>Unused</i>	
4	H2	Geranylgeranylacetone	20218
4	H3	Cholic Acid	20250
4	H4	Lithocholic Acid	20253
4	H5	Glycodeoxycholic Acid (hydrate)	20274
4	H6	Taurochenodeoxycholic Acid (sodium salt)	20275
4	H7	Glycocholic Acid	20276
4	H8	Tauroursodeoxycholic Acid	20277
4	H9	β -Muricholic Acid	20287
4	H10	Murideoxycholic Acid	20290
4	H11	ω -Muricholic Acid	20292
4	H12	<i>Unused</i>	
5	A1	<i>Unused</i>	
5	A2	15(R)-Prostaglandin D ₂	10118
5	A3	24(S),25-epoxy Cholesterol	10131
5	A4	Isopropyl dodec-11-enylfluorophosphonate	11613
5	A5	Prostaglandin D ₂ -1-glycerol ester	12015
5	A6	Prostaglandin E ₁	13010
5	A7	13,14-dihydro-15-keto-tetranor Prostaglandin E ₂	13101
5	A8	9-deoxy-9-methylene-16,16-dimethyl Prostaglandin E ₂	14420
5	A9	11 β -Prostaglandin E ₂	14510
5	A10	4(Z),7(Z),10(Z),13(Z),16(Z)-N-(2-hydroxyethyl)-Docosapentaenamide	14723
5	A11	16,16-dimethyl Prostaglandin E ₂ p-(p-acetamidobenzamido) phenyl ester	14753
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



Page 12 of 35

Plate	Well	Contents	Item Number
5	B1	<i>Unused</i>	
5	B2	16 α -hydroxy Estrone	15208
5	B3	1a,1b-dihomo Prostaglandin F _{2α}	16050
5	B4	11 β -Prostaglandin F _{2α} Ethanolamide	16522
5	B5	9-keto Fluprostenol isopropyl ester	16782
5	B6	5-POHSA	17114
5	B7	Arachidonoyl Chloride	18078
5	B8	Carbaprostacyclin	18210
5	B9	Hyodeoxycholic Acid	20294
5	B10	21-desacetyl Deflazacort	20316
5	B11	Galactosylsphingosine (d18:1)	20338
5	B12	<i>Unused</i>	
5	C1	<i>Unused</i>	
5	C2	Betamethasone	20363
5	C3	N-Oleoyl Alanine	20368
5	C4	N-Oleoyl Proline	20369
5	C5	Deflazacort	20386
5	C6	Estrone 3-sulfate (sodium salt)	20513
5	C7	13(Z)-Docosenoic Acid methyl ester	20568
5	C8	11(Z)-Eicosenoic Acid	20606
5	C9	Lithocholic Acid 3-sulfate (sodium salt)	20676
5	C10	Prednisone	20677
5	C11	Fluticasone Propionate	20703
5	C12	<i>Unused</i>	
5	D1	<i>Unused</i>	
5	D2	13Z,16Z-Docosadienoic Acid	20749
5	D3	Deoxycholic Acid	20756
5	D4	17 α -Estradiol	20776
5	D5	Marinobufagenin	20798
5	D6	Decanoic Acid (sodium salt)	20838
5	D7	Acitretin	20853
5	D8	Flurandrenolide	20933
5	D9	1-Myristoyl-2-hydroxy- <i>sn</i> -glycero-3-PC	20958
5	D10	1-Oleoyl-2-hydroxy- <i>sn</i> -glycero-3-PC	20959
5	D11	1-Oleoyl-2-hydroxy- <i>sn</i> -glycero-3-PG (sodium salt)	20960
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 35

Plate	Well	Contents	Item Number
5	E1	<i>Unused</i>	
5	E2	1-Palmitoyl-2-hydroxy-sn-glycero-3-PG (sodium salt)	20961
5	E3	1-Stearoyl-2-hydroxy-sn-glycero-3-PG (sodium salt)	20964
5	E4	Linoleic Acid Amide	21055
5	E5	Elaidamide	21079
5	E6	Hexadecanamide	21086
5	E7	Stearamide	21087
5	E8	Hecogenin	21372
5	E9	DCP-LA	21375
5	E10	22(S)-hydroxy Cholesterol	21399
5	E11	10(S)-PAHSA	21419
5	E12	<i>Unused</i>	
5	F1	<i>Unused</i>	
5	F2	Brassinolide	21594
5	F3	Glycoursodeoxycholic Acid	21698
5	F4	Glycolithocholic Acid	21723
5	F5	Cholesteryl Heptadecanoate	21750
5	F6	Fucosterol	21858
5	F7	3 β -hydroxy-5-Cholestenoic Acid	21859
5	F8	Taurohyodeoxycholic Acid (sodium salt)	21956
5	F9	β -Estradiol 17-valerate	22221
5	F10	Brassicasterol	22296
5	F11	Myristelaidic Acid	22361
5	F12	<i>Unused</i>	
5	G1	<i>Unused</i>	
5	G2	10(Z)-Pentadecenoic Acid	22362
5	G3	10(E)-Pentadecenoic Acid	22467
5	G4	<i>trans</i> -10-Heptadecenoic Acid	22468
5	G5	Ricinelaiddic Acid	22470
5	G6	Aramchol	22478
5	G7	Sphinganine-1-phosphate (d18:0)	22500
5	G8	N,N-dimethyl Sphingosine (d17:1)	22503
5	G9	N,N-dimethyl Sphinganine (d18:0)	22504
5	G10	C4 Ceramide (d18:1/4:0)	22528
5	G11	C10 Ceramide (d18:1/10:0)	22529
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	C12 Ceramide (d18:1/12:0)	22530
5	H3	C14 Ceramide (d18:1/14:0)	22531
5	H4	C16 Ceramide-1-phosphate (d18:1/16:0) (ammonium salt)	22542
5	H5	8(Z)-Eicosenoic Acid	22579
5	H6	13(E)-Docosenoic Acid	22580
5	H7	Heneicosanoic Acid	22593
5	H8	12(Z),15(Z)-Heneicosadienoic Acid	22594
5	H9	Cholesteryl Arachidonate	22595
5	H10	Cholesterol α -Linolenate	22596
5	H11	Cholesteryl Linoleate	22597
5	H12	<i>Unused</i>	
6	A1	<i>Unused</i>	
6	A2	Glycohyodeoxycholic Acid	22643
6	A3	Anacardic Acid diene	22663
6	A4	3-hydroxy Docosanoic Acid	22666
6	A5	3-hydroxy Lignoceric Acid	22667
6	A6	Taurohyocholic Acid (sodium salt)	22669
6	A7	Glycohyocholic Acid	22670
6	A8	2-hydroxy Decanoic Acid	22678
6	A9	2-hydroxy Palmitic Acid	22679
6	A10	2-hydroxy Lignoceric Acid	22680
6	A11	C18 Phytoceramide (t18:0/18:0)	22686
6	A12	<i>Unused</i>	
6	B1	<i>Unused</i>	
6	B2	C18 (($\pm$)-2'-hydroxy) Ceramide (d18:1/18:0)	22687
6	B3	3-hydroxy Tridecanoic Acid	22689
6	B4	3-hydroxy Stearic Acid	22692
6	B5	Cholesteryl Docosapentaenoate	22718
6	B6	Cholesteryl Eicosapentaenoate	22719
6	B7	C2 Phytoceramide (t18:0/2:0)	22823
6	B8	3-keto Sphinganine (d8:0) (hydrochloride)	22824
6	B9	C24 Phytoceramide (t18:0/24:0)	22827
6	B10	Oleoyl Ethanolamide Phosphate	22886
6	B11	11-deoxy Corticosterone	22916
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	9(E),11(E),13(E)-Octadecatrienoic Acid	22976
6	C3	Estramustine Phosphate (sodium salt)	23126
6	C4	C22 Glucosylceramide (d18:1/22:0)	23210
6	C5	1-β-D-Glucosylsphingosine (d18:1)	23211
6	C6	1-β-D-Glucosylsphingadienine (d18:2 (4E,8E))	23213
6	C7	Equilenin	23322
6	C8	Drospirenone	23347
6	C9	20(S)-Ginsenoside Rh ₂	23390
6	C10	4(Z),7(Z)-Decadienoic Acid	23516
6	C11	Desogestrel	23651
6	C12	<i>Unused</i>	
6	D1	<i>Unused</i>	
6	D2	Medroxyprogesterone 17-acetate	23664
6	D3	3-Formylsalicylic Acid	23368
6	D4	6-Aminocaproic Acid	23724
6	D5	Halobetasol Propionate	23781
6	D6	Docosanol	23793
6	D7	Cortisone Acetate	23798
6	D8	Fluocinonide	23803
6	D9	Estropipate	23806
6	D10	Diflorasone Diacetate	23808
6	D11	Mestranol	23822
6	D12	<i>Unused</i>	
6	E1	<i>Unused</i>	
6	E2	Flunisolide	23842
6	E3	CAY10726	23859
6	E4	Alclometasone Dipropionate	23901
6	E5	Amcinonide	23903
6	E6	Desoximetasone	23945
6	E7	Azelaic Acid	23977
6	E8	Hydroxyprogesterone Caproate	23978
6	E9	Flumethasone	24184
6	E10	Sulfatides (bovine) (sodium salt)	24323
6	E11	C8 Galactosylceramide (d18:1/8:0)	24346
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



Page 16 of 35

Plate	Well	Contents	Item Number
6	F1	<i>Unused</i>	
6	F2	C12 Galactosylceramide (d18:1/12:0)	24347
6	F3	C16 Lactosylceramide (d18:1/16:0)	24352
6	F4	C8 Phytoceramide (t18:0/8:0)	24357
6	F5	DL-erythro/threo Sphinganine (d18:0)	24367
6	F6	C6 dihydro Ceramide (d18:0/6:0)	24368
6	F7	L-erythro Sphingosine (d18:1)	24372
6	F8	L-erythro Sphinganine (d18:0)	24374
6	F9	D-threo Sphinganine (d18:0)	24375
6	F10	DL-erythro/threo Sphinganine (d16:0)	24376
6	F11	DL-erythro Sphinganine (d20:0)	24378
6	F12	<i>Unused</i>	
6	G1	<i>Unused</i>	
6	G2	C18 dihydro Ceramide (d18:0/18:0)	24379
6	G3	3-keto Sphinganine (d12:0) (hydrochloride)	24384
6	G4	N,N-dihexyl Sphingosine (d18:1/6:0/6:0)	24385
6	G5	C2 L-threo Ceramide (d18:1/2:0)	24386
6	G6	C6 L-erythro Ceramide (d18:1/6:0)	24388
6	G7	C6 L-threo Ceramide (d18:1/6:0)	24390
6	G8	C18:1 Ceramide (d18:1/18:1(9Z))	24397
6	G9	C32 Ceramide (d18:1/32:0)	24424
6	G10	C2 Ceramide (d14:1/2:0)	24425
6	G11	C12 ((±)-2'-hydroxy) dihydro Ceramide (d18:0/12:0)	24435
6	G12	<i>Unused</i>	
6	H1	<i>Unused</i>	
6	H2	C18 ((±)-2'-hydroxy) dihydro Ceramide (d18:0/18:0)	24437
6	H3	L-threo Lysosphingomyelin (d18:1)	24454
6	H4	Lysodihydrosphingomyelin (d18:0)	24463
6	H5	C2 Galactosylceramide (d18:1/2:0)	24465
6	H6	C15 Galactosylceramide (d18:1/15:0)	24466
6	H7	C6 Glucosylceramide (d18:2/6:0)	24474
6	H8	1-Deoxysphingosine (m18:1(14Z))	24515
6	H9	2-fluoro Palmitic Acid methyl ester	24526
6	H10	9(Z),11(Z)-Conjugated Linoleic Acid	24582
6	H11	2-hydroxy Lauric Acid	24586
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



Page 17 of 35

Plate	Well	Contents	Item Number
7	A1	<i>Unused</i>	
7	A2	24-dehydro Cholesterol	14943
7	A3	Lanosterol	19521
7	A4	2-hydroxy Stearic Acid	24590
7	A5	2-hydroxy Arachidic Acid	24592
7	A6	3-hydroxy Octanoic Acid	24609
7	A7	3-hydroxy Decanoic Acid	24613
7	A8	N-Glycine Galactosylsphingosine (d18:1)	24620
7	A9	3-hydroxy Undecanoic Acid	24640
7	A10	3-hydroxy Lauric Acid	24642
7	A11	3-hydroxy Myristic Acid methyl ester	24645
7	A12	<i>Unused</i>	
7	B1	<i>Unused</i>	
7	B2	20-hydroxy Arachidic Acid	24652
7	B3	22-hydroxy Docosanoic Acid	24655
7	B4	1-Deoxysphingosine (m18:1(4E))	24687
7	B5	12-methyl Tridecanoic Acid methyl ester	24810
7	B6	13-methyl Myristic Acid	24811
7	B7	14-methyl Pentadecanoic Acid methyl ester	24813
7	B8	15-methyl Palmitic Acid	24814
7	B9	12-methyl Myristic Acid	24817
7	B10	13-methyl Pentadecanoic Acid methyl ester	24819
7	B11	14-methyl Palmitic Acid	24820
7	B12	<i>Unused</i>	
7	C1	<i>Unused</i>	
7	C2	DL-2,6-dimethyl Heptanoic Acid	24822
7	C3	10-methyl Palmitic Acid methyl ester	24823
7	C4	cis-9,10-Methyleneoctadecanoic Acid	24824
7	C5	Lactosylsphingosine (d18:1) (synthetic)	24867
7	C6	Globotetraosylceramides (porcine RBC)	24881
7	C7	Medroxyprogesterone	24908
7	C8	Ginkgolide A	24970
7	C9	17-phenyl trinor Prostaglandin F _{2α} dimethyl amide	25241
7	C10	17-phenyl trinor Prostaglandin F _{2α} isopropyl amide	25242
7	C11	Photoclick Sphingosine	25364
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	1a-monohomo-17-phenyl trinor Prostaglandin F _{2α}	25392
7	D3	7α-hydroxy-4-Cholesten-3-one	25442
7	D4	17-phenyl trinor Prostaglandin F _{2α} glycinate	25470
7	D5	C6 Urea Ceramide	25491
7	D6	C6 1-Deoxyceramide (m18:1(14Z)/6:0)	25493
7	D7	C6 dihydro 1-Deoxyceramide (m18:0/6:0)	25494
7	D8	5α,6β-Dihydroxycholesterol	25538
7	D9	5α,6α-epoxy Cholesterol	25539
7	D10	27-hydroxy Cholestenone	25550
7	D11	1-Myristoyl-2-hydroxy-sn-glycero-3-PG (sodium salt)	25587
7	D12	<i>Unused</i>	
7	E1	<i>Unused</i>	
7	E2	1-Myristoyl-2-hydroxy-sn-glycero-3-PE	25591
7	E3	1-Stearoyl-2-hydroxy-sn-glycero-3-PE	25593
7	E4	1-Oleoyl-2-hydroxy-sn-glycero-3-PE	25596
7	E5	C24 dihydro Ceramide (d18:0/24:0)	25602
7	E6	5β,6β-epoxy Cholesterol	25603
7	E7	Betamethasone 17-valerate	25640
7	E8	1-Palmitoleoyl glycerol	25697
7	E9	Cholesteryl Hemisuccinate	25698
7	E10	6α-hydroxy Cholesterol	25715
7	E11	1-O-Hexadecyl-sn-glycerol	25723
7	E12	<i>Unused</i>	
7	F1	<i>Unused</i>	
7	F2	Dihexadecyl Phosphate	25727
7	F3	Glucosylceramide (plant)	25850
7	F4	12,15-epoxy-13-methyl-12,14-Eicosadienoic Acid	25862
7	F5	N-[1-(2,3-Dioleoyloxy)propyl]-N,N,N-trimethylammonium (chloride)	25926
7	F6	Cho-Arg (trifluoroacetate salt)	25943
7	F7	2H-Cho-Arg (trifluoroacetate salt)	25944
7	F8	Dios-Arg (trifluoroacetate salt)	25945
7	F9	8-dehydro Cholesterol	25970
7	F10	7-keto-25-hydroxy Cholesterol	25973
7	F11	Hecogenin Acetate	25977
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	Heptadecanoyl-Coenzyme A (lithium salt hydrate)	25999
7	G3	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PE	26011
7	G4	9(Z),11(E),13(Z)-Octadecatrienoic Acid	26057
7	G5	Sitostanol	26094
7	G6	Trihydroxycholestanoic Acid	26098
7	G7	2-chloro Palmitic Acid	26107
7	G8	Cholic Acid anilide	26144
7	G9	13(Z)-Eicosenoic Acid	26169
7	G10	Zymostenol	26209
7	G11	Cholesteryl Myristate	26439
7	G12	<i>Unused</i>	
7	H1	<i>Unused</i>	
7	H2	Cholesteryl Arachidate	26465
7	H3	Cholesteryl Behenate	26466
7	H4	Cholesteryl 11,14-Eicosadienoate	26467
7	H5	Cholesteryl Erucate	26468
7	H6	Cholesteryl Dihomo- γ -Linolenate	26469
7	H7	Cholesteryl Lignocerate	26470
7	H8	Cholesteryl Oleate	26472
7	H9	Cholesteryl Palmitate	26473
7	H10	Cholesteryl Palmitoleate	26474
7	H11	Cholesteryl Stearate	26475
7	H12	<i>Unused</i>	
8	A1	<i>Unused</i>	
8	A2	Carbocyclic Thromboxane A ₂	19010
8	A3	3-keto Sphinganine (d18:0) (hydrochloride)	24380
8	A4	2-hydroxy Docosanoic Acid	24594
8	A5	Tetrahydro-11-deoxy Cortisol	26501
8	A6	Campestanol	26538
8	A7	Oleoyl-L-carnitine	26557
8	A8	Myristoyl-L-carnitine (chloride)	26559
8	A9	Linoleoyl-L-carnitine (chloride)	26560
8	A10	DMHAPC-Chol	26582
8	A11	DMPAC-Chol	26583
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	HAPC-Chol	26584
8	B3	MHAPC-Chol	26585
8	B4	OH-Chol	26586
8	B5	tetranor-Prostaglandin E ₁	26702
8	B6	Cholesterol β-D-Glucoside	26713
8	B7	11-Hexadecenoic Acid	26730
8	B8	Sterculic Acid	26735
8	B9	5α-Cholestane	26763
8	B10	Coprostanol	26764
8	B11	17-methyl Stearic Acid	26766
8	B12	<i>Unused</i>	
8	C1	<i>Unused</i>	
8	C2	1-Oleoyl Glycerol	26793
8	C3	1-Stearoyl- <i>rac</i> -glycerol	26827
8	C4	Tricosanoic Acid	26867
8	C5	Monodecanoin	26888
8	C6	Sphingosine (d12:1)	26893
8	C7	1-Octanoyl- <i>rac</i> -glycerol	26945
8	C8	1-Arachidoyl- <i>rac</i> -glycerol	26947
8	C9	1-Pentadecanoyl- <i>rac</i> -glycerol	26959
8	C10	1-Tridecanoyl- <i>rac</i> -glycerol	26961
8	C11	1-Palmitoyl- <i>rac</i> -glycerol	26993
8	C12	<i>Unused</i>	
8	D1	<i>Unused</i>	
8	D2	C22:1 1-Deoxyceramide (m18:1/22:1)	27015
8	D3	C22 dihydro 1-Deoxyceramide (m18:0/22:0)	27018
8	D4	Sphingosine (d18:1(14Z))	27019
8	D5	Sphingosine-1-phosphate (d18:1(14Z))	27020
8	D6	Squalene	27058
8	D7	16α-hydroxy Prednisolone	27169
8	D8	Lactosylceramides (bovine buttermilk)	27197
8	D9	Esterified Steryl Glucosides	27204
8	D10	Steryl Glucosides	27205
8	D11	Altrenogest	27218
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

Plate	Well	Contents	Item Number
8	E1	<i>Unused</i>	
8	E2	2-Stearoyl- <i>rac</i> -glycerol	27282
8	E3	Dihexanoin	27293
8	E4	(R)- α -Lipoic Acid	27299
8	E5	Ginsenoside Rh ₁	27318
8	E6	<i>cis</i> -8-Octadecenoic Acid	27447
8	E7	C22 1-Deoxyceramide (m18:1(14Z)/22:0)	27561
8	E8	C24:1 dihydro 1-Deoxyceramide (m18:0/24:1)	27562
8	E9	5 α -hydroxy Laxogenin	27665
8	E10	16 α -Hydroxyetiocholanolone	27683
8	E11	Ganoderenic Acid A	27685
8	E12	<i>Unused</i>	
8	F1	<i>Unused</i>	
8	F2	Butyryl-Coenzyme A (sodium salt)	27865
8	F3	Hexanoyl-Coenzyme A (sodium salt)	27866
8	F4	Octanoyl-Coenzyme A (sodium salt)	27868
8	F5	Estrone 3-methyl ether	27978
8	F6	Telocinobufagin	27981
8	F7	17 β -hydroxy Exemestane	27989
8	F8	1-Octacosanol	28010
8	F9	2-O-Acetyl-20-hydroxyecdysone	28063
8	F10	Solasonine	28111
8	F11	1-Lauroyl- <i>rac</i> -glycerol	28170
8	F12	<i>Unused</i>	
8	G1	<i>Unused</i>	
8	G2	1-Myristoyl- <i>rac</i> -glycerol	28173
8	G3	1-Heptadecanoyl- <i>rac</i> -glycerol	28180
8	G4	15(S)-acetate Prostaglandin A ₂ methyl ester	28195
8	G5	Glyco-Obeticholic Acid	28242
8	G6	Tauro-Obeticholic Acid	28243
8	G7	Triamcinolone	28275
8	G8	3 α ,5 α -THDOC	28364
8	G9	3-Oxochoolic Acid	28403
8	G10	12-hydroxy Lauric Acid	28494
8	G11	C16 Ceramide (d18:1(14Z)/16:0)	28601
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



Page 22 of 35

Plate	Well	Contents	Item Number
8	H1	<i>Unused</i>	
8	H2	C22:1 Ceramide (d18:1(14Z)/22:1(13Z))	28603
8	H3	C24 Ceramide (d18:1(14Z)/24:0)	28604
8	H4	Decanoyl-Coenzyme A (hydrate)	28607
8	H5	Lauroyl-Coenzyme A (potassium salt hydrate)	28610
8	H6	9,10-Dihydroxystearic Acid	28612
8	H7	17R(18S)-EpETE	28630
8	H8	3-hydroxy Heptadecanoic Acid	28735
8	H9	1-Undecanoyl- <i>rac</i> -glycerol	28763
8	H10	Tauro- ω -muricholic Acid (sodium salt)	28842
8	H11	N-Oleoyl Glutamine	28922
8	H12	<i>Unused</i>	
9	A1	<i>Unused</i>	
9	A2	C24 (2'(R)-hydroxy) Phytoceramide (t18:0/24:0)	28933
9	A3	C24 (2'(R)-hydroxy) dihydro Ceramide (d18:0/24:0)	28935
9	A4	Ginsenoside Rb ₃	29005
9	A5	20(S)-Protopanaxatriol	29073
9	A6	Dehydrocholic Acid	29176
9	A7	17 α -hydroxy Pregnenolone	29181
9	A8	Hydrocortisone Acetate	29390
9	A9	Fludrocortisone	29391
9	A10	Chlormadinone Acetate	29392
9	A11	pNPS-DHA	29507
9	A12	<i>Unused</i>	
9	B1	<i>Unused</i>	
9	B2	14-dehydro Zymostenol	29534
9	B3	3 β -OH-7-Oxocholeic Acid	29541
9	B4	Alloisolithocholic Acid	29542
9	B5	Choleic Acid	29543
9	B6	Dehydrolithocholic Acid	29544
9	B7	Isolithocholic Acid	29545
9	B8	7-keto Lithocholic Acid	29546
9	B9	12-keto Lithocholic Acid	29547
9	B10	Hentriacontane	29776
9	B11	Betamethasone 21-phosphate (sodium salt hydrate)	29778
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



Page 23 of 35

Plate	Well	Contents	Item Number
9	C1	<i>Unused</i>	
9	C2	Ethynodiol Diacetate	29854
9	C3	Fluticasone Furoate	29878
9	C4	9-Oxononanoic Acid	29882
9	C5	Ursocholic Acid	29883
9	C6	Isodeoxycholic Acid	29890
9	C7	Glycodehydrocholic Acid	29892
9	C8	Ginsenoside CK	29926
9	C9	C6 1-Deoxyceramide (m18:1/6:0)	30027
9	C10	C12 1-Deoxyceramide (m18:1/12:0)	30028
9	C11	C16 1-Deoxyceramide (m18:1/16:0)	30029
9	C12	<i>Unused</i>	
9	D1	<i>Unused</i>	
9	D2	C24:1 1-Deoxyceramide (m18:1/24:1(15Z))	30030
9	D3	Myristoyl Glycine	30107
9	D4	Tigogenin	30137
9	D5	Δ^5 -Avenasterol	30154
9	D6	9(E),11(E)-9-nitro Conjugated Linoleic Acid	30160
9	D7	Oleyl Alcohol	30385
9	D8	Allocholic Acid	30415
9	D9	14-Anhydrodigitoxigenin	30439
9	D10	10-hydroxy Decanoic Acid	30458
9	D11	Elaidyl Alcohol	30459
9	D12	<i>Unused</i>	
9	E1	<i>Unused</i>	
9	E2	12-Bromododecanoic Acid	30465
9	E3	Arenobufagin	30695
9	E4	Cortisone	30763
9	E5	Nordeoxycholic Acid	30837
9	E6	Deoxycorticosterone Acetate	30856
9	E7	Ergosterol Peroxide	30860
9	E8	21-Deoxycortisol	30873
9	E9	8(E),10(E),12(Z)-Octadecatrienoic Acid	30885
9	E10	Nomegestrol	30893
9	E11	ATRA-BA Hybrid	30982
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

Plate	Well	Contents	Item Number
9	F1	<i>Unused</i>	
9	F2	11(Z),14(Z)-Eicosadienoic Acid Ethanolamide	31024
9	F3	Estradiol Cypionate	31105
9	F4	5 α -Dihydroprogesterone	31505
9	F5	4-Hydroxyestradiol	31508
9	F6	Mometasone	31702
9	F7	Octacosanoic Acid	31735
9	F8	4-hydroxy Hexenal	32060
9	F9	4-hydroxy Nonenal	32100
9	F10	4-hydroxy Nonenal Mercapturic Acid	32110
9	F11	($\pm$)9-HODE cholesteryl ester	38401
9	F12	<i>Unused</i>	
9	G1	<i>Unused</i>	
9	G2	($\pm$)5(6)-DiHET	51211
9	G3	($\pm$)5(6)-DiHET lactone	51212
9	G4	PAF C-16	60900
9	G5	1-O-hexadecyl-2-Dihomo- γ -Linolenoyl- <i>sn</i> -glycero-3-PC	60901
9	G6	2-O-methyl PAF C-16	60902
9	G7	1-O-hexadecyl-2-Docosahexaenoyl- <i>sn</i> -glycero-3-PC	60903
9	G8	1-O-hexadecyl-2-Arachidonoyl- <i>sn</i> -glycero-3-PC	60904
9	G9	Hexanolamino PAF C-16	60905
9	G10	Lyso-PAF C-16	60906
9	G11	1-O-hexadecyl-2-Eicosapentaenoyl- <i>sn</i> -glycero-3-PC	60907
9	G12	<i>Unused</i>	
9	H1	<i>Unused</i>	
9	H2	Methylcarbamoyl PAF C-16	60908
9	H3	Pyrrolidino PAF C-16	60909
9	H4	PAF C-18	60910
9	H5	2-O-methyl PAF C-18 (hydrate)	60912
9	H6	Lyso-PAF C-18	60916
9	H7	Hexadecyl Acetyl Glycerol	60920
9	H8	Azelaoyl PAF	60924
9	H9	2-O-ethyl PAF C-16	60925
9	H10	Butanoyl PAF	60928
9	H11	Butenoyl PAF	60929
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 35

Plate	Well	Contents	Item Number
10	A1	<i>Unused</i>	
10	A2	Hexadecyl Methyl Glycerol	60930
10	A3	2-thio PAF	60945
10	A4	Arachidonyl Trifluoromethyl Ketone	62120
10	A5	1-Arachidonoyl Glycerol	62150
10	A6	2-Arachidonoyl Glycerol	62160
10	A7	2-Arachidonoyl Glycerol ether	62165
10	A8	Arachidonoyl Serinol	62170
10	A9	Palmitoyl Serinol	62175
10	A10	1-Oleoyl-2-hydroxy-sn-glycero-3-PA (sodium salt)	62215
10	A11	1,2-bis(heptanoylthio) Glycerophosphocholine	62235
10	A12	<i>Unused</i>	
10	B1	<i>Unused</i>	
10	B2	2-Linoleoyl Glycerol	62260
10	B3	C2 Ceramide (d18:1/2:0)	62510
10	B4	C6 Ceramide (d18:1/6:0)	62525
10	B5	C24:1 Ceramide (d18:1/24:1(15Z))	62530
10	B6	C24 Ceramide (d18:1/24:0)	62535
10	B7	C8 Ceramine (d18:1/8:0)	62545
10	B8	C8 Ceramide-1-phosphate (d18:1/8:0)	62547
10	B9	Sphingosine-1-phosphate (d18:1)	62570
10	B10	N,N-Dimethylsphingosine (d18:1)	62575
10	B11	Oleoyl Trifluoromethyl Ketone	62640
10	B12	<i>Unused</i>	
10	C1	<i>Unused</i>	
10	C2	Thioetheramide-PC	62750
10	C3	7-hydroxycoumarinyl Arachidonate	62910
10	C4	PAz-PC	62924
10	C5	KDdiA-PC	62935
10	C6	KOdiA-PC	62945
10	C7	DMAPP (ammonium salt)	63180
10	C8	Farnesyl Pyrophosphate (ammonium salt)	63250
10	C9	S-Farnesyl Thioacetic Acid	63260
10	C10	N-acetyl-S-farnesyl-L-Cysteine	63270
10	C11	Geranyl Pyrophosphate (triammonium salt)	63320
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



Page 26 of 35

Plate	Well	Contents	Item Number
10	D1	<i>Unused</i>	
10	D2	Geranylgeranyl Pyrophosphate (ammonium salt)	63330
10	D3	α -hydroxy Farnesyl Phosphonic Acid	63420
10	D4	Lycopene	70945
10	D5	<i>cis</i> -Parinaric Acid	71430
10	D6	($\pm$)-Jasmonic Acid	88300
10	D7	12-oxo Phytodienoic Acid	88520
10	D8	Traumatic Acid	88820
10	D9	n-Triacontanol	88840
10	D10	22(R)-hydroxy Cholesterol	89355
10	D11	Arachidonic Acid	90010
10	D12	<i>Unused</i>	
10	E1	<i>Unused</i>	
10	E2	ω -3 Arachidonic Acid	90011
10	E3	5,6-dehydro Arachidonic Acid	90020
10	E4	Arachidonoyl Ethanolamide	90050
10	E5	Arachidonoyl Glycine	90051
10	E6	Arvanil	90052
10	E7	Arachidonoyl-2'-Fluoroethylamide	90054
10	E8	($\pm$)-2-Methyl Arachidonoyl-2'-Fluoroethylamide	90055
10	E9	N-Arachidonoyl Dopamine	90057
10	E10	Arachidonoyl <i>m</i> -Nitroaniline	90059
10	E11	AM404	90060
10	E12	<i>Unused</i>	
10	F1	<i>Unused</i>	
10	F2	N-Arachidonoyl-L-Alanine	90065
10	F3	N-Arachidonoyl- γ -Aminobutyric Acid	90067
10	F4	R-1 Methanandamide	90070
10	F5	S-1 Methanandamide	90072
10	F6	R-2 Methanandamide	90074
10	F7	S-2 Methanandamide	90076
10	F8	Eicosapentaenoic Acid	90110
10	F9	5,8,11,14-Eicosatetraenoic Acid	90120
10	F10	9(Z),11(E)-Conjugated Linoleic Acid	90140
10	F11	10(E),12(Z)-Conjugated Linoleic Acid	90145
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



Page 27 of 35

Plate	Well	Contents	Item Number
10	G1	<i>Unused</i>	
10	G2	Linoleic Acid	90150
10	G3	Linoleoyl Ethanolamide	90155
10	G4	Linoelaidic Acid	90160
10	G5	Docosapentaenoic Acid	90165
10	G6	Docosatrienoic Acid	90170
10	G7	13(Z)-Docosenoic Acid	90175
10	G8	5(Z),8(Z),11(Z)-Eicosatrienoic Acid	90190
10	G9	11(Z),14(Z),17(Z)-Eicosatrienoic Acid	90192
10	G10	5(Z),8(Z),11(Z)-Eicosatrienoic Acid Ethanolamide	90195
10	G11	5,8,11-Eicosatriynoic Acid	90200
10	G12	<i>Unused</i>	
10	H1	<i>Unused</i>	
10	H2	α -Linolenic Acid	90210
10	H3	γ -Linolenic Acid	90220
10	H4	Dihomo- γ -Linolenic Acid	90230
10	H5	Dihomo- γ -Linolenoyl Ethanolamide	90235
10	H6	Elaidic Acid	90250
10	H7	Oleic Acid	90260
10	H8	Olvanil	90262
10	H9	Oleoyl Ethanolamide	90265
10	H10	N-Oleoyl Glycine	90269
10	H11	17-Octadecynoic Acid	90270
10	H12	<i>Unused</i>	
11	A1	<i>Unused</i>	
11	A2	MEDICA 16	90290
11	A3	Adrenic Acid	90300
11	A4	Docosahexaenoic Acid	90310
11	A5	Stearidonic Acid	90320
11	A6	11(Z),14(Z)-Eicosadienoic Acid	90330
11	A7	Heptadecanoyl Ethanolamide	90342
11	A8	Palmitoyl Ethanolamide	90350
11	A9	R-Palmitoyl-(1-methyl) Ethanolamide	90353
11	A10	R-Palmitoyl-(2-methyl) Ethanolamide	90357
11	A11	Phytanic Acid	90360
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



Page 28 of 35

Plate	Well	Contents	Item Number
11	B1	<i>Unused</i>	
11	B2	9(E),11(E)-Conjugated Linoleic Acid	90370
11	B3	Oleamide	90375
11	B4	13-Docosenamide	90377
11	B5	2-fluoro Palmitic Acid	90380
11	B6	Docosatetraenoyl Ethanolamide	90385
11	B7	2-hydroxy Myristic Acid	90390
11	B8	9,12-Octadecadiynoic Acid	90400
11	B9	3-Thiatetradecanoic Acid	90500
11	B10	O-11	90510
11	B11	O-Arachidonoyl Ethanolamine (hydrochloride)	91050
11	B12	<i>Unused</i>	
11	C1	<i>Unused</i>	
11	C2	Arachidonoyl Cyclopropylamide	91053
11	C3	Arachidonoyl 2'-Chloroethylamide	91054
11	C4	Palmitoyl N-Isopropylamide	91354
11	C5	Petromyzonol	98250
11	C6	Δ^{14} -Triamcinolone acetonide	9000087
11	C7	15-keto Prostaglandin A ₁	9000142
11	C8	Fluorometholone	9000157
11	C9	8-methyl Nonanoic Acid	9000310
11	C10	7(Z)-Tricosene	9000313
11	C11	7(Z),11(Z)-Nonacosadiene	9000314
11	C12	<i>Unused</i>	
11	D1	<i>Unused</i>	
11	D2	Arachidoyl glycine	9000325
11	D3	Linoleoyl Glycine	9000326
11	D4	Stearidonoyl Glycine	9000327
11	D5	Docosahexaenoyl Glycine	9000328
11	D6	Methylcarbaryl PAF C-8	9000332
11	D7	Docosanoic Acid	9000338
11	D8	Arachidic Acid	9000339
11	D9	Oxy-16	9000395
11	D10	13,14-dihydro-16,16-difluoro Prostaglandin E ₁	9000405
11	D11	7(Z)-Pentacosene	9000530
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 35

Plate	Well	Contents	Item Number
11	E1	<i>Unused</i>	
11	E2	7(Z),11(Z)-Pentacosadiene	9000531
11	E3	Oleoyl Serotonin	9000629
11	E4	Palmitoyl Serotonin	9000630
11	E5	Docosahexaenoyl Serotonin	9000639
11	E6	Eicosapentaenoyl Serotonin	9000640
11	E7	17-phenyl trinor Prostaglandin F _{2α} cyclohexylamide	9000686
11	E8	C10 AV Ceramide (d18:1/10:0)	9000753
11	E9	16-hydroxy Hexadecanoic Acid	9000789
11	E10	Tafluprost ethyl amide	9000843
11	E11	Arachidonoyl-L-carnitine (chloride)	9000939
11	E12	<i>Unused</i>	
11	F1	<i>Unused</i>	
11	F2	(R)-(+)-Eicosapentaenyl-1'-Hydroxy-2'-Propylamide	9001225
11	F3	(R)-(-)-Eicosapentaenyl-2'-Hydroxy-1'-Propylamide	9001226
11	F4	(S)-(-)-Eicosapentaenyl-1'-Hydroxy-2'-Propylamide	9001227
11	F5	(S)-(+)-Eicosapentaenyl-2'-Hydroxy-1'-Propylamide	9001228
11	F6	(R)-(+)-Docosahexaenyl-1'-Hydroxy-2'-Propylamide	9001229
11	F7	(R)-(-)-Docosahexaenyl-2'-Hydroxy-1'-Propylamide	9001230
11	F8	(S)-(+)-Docosahexaenyl-2'-Hydroxy-1'-Propylamide	9001231
11	F9	(S)-(-)-Docosahexaenyl-1'-Hydroxy-2'-Propylamide	9001232
11	F10	(R)-(+)-Linoleyl-1'-Hydroxy-2'-Propylamide	9001233
11	F11	(R)-(-)-Linoleyl-2'-Hydroxy-1'-Propylamide	9001234
11	F12	<i>Unused</i>	
11	G1	<i>Unused</i>	
11	G2	(S)-(+)-Linoleyl-2'-Hydroxy-1'-Propylamide	9001235
11	G3	(S)-(-)-Linoleyl-1'-Hydroxy-2'-Propylamide	9001236
11	G4	Resibufogenin	9001348
11	G5	N-Docosahexaenoyl Dopamine	9001394
11	G6	Pentadecanoyl Ethanolamide	9001740
11	G7	Myristoyl Ethanolamide	9001742
11	G8	Tricosanoyl Ethanolamide	9001743
11	G9	Lignoceroyl Ethanolamide	9001744
11	G10	Docosaenoyl Ethanolamide	9001745
11	G11	Nervonoyl Ethanolamide	9001746
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



Page 30 of 35

Plate	Well	Contents	Item Number
11	H1	<i>Unused</i>	
11	H2	γ -Linolenoyl Ethanolamide	9001747
11	H3	Palmitelaidic Acid	9001798
11	H4	Linoleoyl Ethanolamide Phosphate	9001837
11	H5	<i>cis</i> -6-Hexadecenoic Acid	9001845
11	H6	Propionyl-L-carnitine (chloride)	9001873
11	H7	Arachidonic Acid sulfonate (sodium salt)	9001886
11	H8	Hexadecanal	9001996
11	H9	4(Z),7(Z),10(Z),13(Z)-Hexadecatetraenoic Acid	9002202
11	H10	N-Benzylpalmitamide	9002235
11	H11	CAY10695	9002312
11	H12	<i>Unused</i>	
12	A1	<i>Unused</i>	
12	A2	CAY10696	9002313
12	A3	N-(8Z-Heptadecen-1-yl)-O-(3-pyridylmethyl) carbamate	9002365
12	A4	Prostaglandin D ₂ Dopamine	9002504
12	A5	Cholic Acid 7-sulfate	9002532
12	A6	Abiraterone	9002768
12	A7	HE-3235	9002769
12	A8	9(Z)-Pentacosene	9002807
12	A9	1-Palmitoyl-2-formyllyl PC	9002943
12	A10	N-(1-(3,4-Dihydroxyphenyl)propan-2-yl)oleamide	9002954
12	A11	<i>Unused</i>	
12	A12	<i>Unused</i>	
12	B1	<i>Unused</i>	
12	B2	Cholesterol	9003100
12	B3	Cholestanol	9003101
12	B4	Lathosterol	9003102
12	B5	Glycine- β -muricholic Acid	9003230
12	B6	Norhyodeoxycholic Acid	9003440
12	B7	Aldosterone 21-sulfate (sodium salt)	9003450
12	B8	Palmitoylcholine (chloride)	9003456
12	B9	C22 Galactosylceramide (d18:1/22:0)	9003458
12	B10	Fluprostenol serinol amide	10004236
12	B11	17-phenyl trinor Prostaglandin F _{2α} serinol amide	10004237
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	<i>trans</i> -4,5-epoxy-2(E)-Decenal	10004257
12	C3	R-1 Methanandamide Phosphate	10004281
12	C4	14,15-EE-5(Z)-E	10004946
12	C5	(+)-5- <i>trans</i> Cloprostenol	10004970
12	C6	5(S),6(R)-7-trihydroxymethyl Heptanoate	10005032
12	C7	AMC Arachidonoyl Amide	10005098
12	C8	9(E),12(Z)-Octadecadienoic Acid	10005146
12	C9	9(Z),12(Z),15(Z),18(Z),21(Z)-Tetracosapentaenoic Acid	10005157
12	C10	6(Z),9(Z),12(Z),15(Z),18(Z),21(Z)-Tetracosahexaenoic Acid	10005165
12	C11	DAUDA	10005188
12	C12	<i>Unused</i>	
12	D1	<i>Unused</i>	
12	D2	AM1172	10005223
12	D3	Tafluprost (free acid)	10005439
12	D4	Tafluprost	10005440
12	D5	N-Arachidonoyl-L-Serine	10005455
12	D6	Oleoyl Ethyl Amide	10005459
12	D7	N-Arachidonoyl Taurine	10005537
12	D8	N-Oleoyl Taurine	10005609
12	D9	N-Palmitoyl Taurine	10005611
12	D10	CAY10448	10005633
12	D11	DL- α -Lipoic Acid	10005728
12	D12	<i>Unused</i>	
12	E1	<i>Unused</i>	
12	E2	Arachidoyl Ethanolamide	10005765
12	E3	Docosanoyl Ethanolamide	10005823
12	E4	17 β -Estradiol	10006315
12	E5	Mifepristone	10006317
12	E6	Levonorgestrel	10006318
12	E7	Norgestrel	10006319
12	E8	Quinestrol	10006320
12	E9	Tibolone	10006321
12	E10	Estriol	10006484
12	E11	Estrone	10006485
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 35

Plate	Well	Contents	Item Number
12	F1	<i>Unused</i>	
12	F2	Ethynyl Estradiol	10006486
12	F3	Estradiol Benzoate	10006487
12	F4	Gestrinone	10006488
12	F5	Thromboxane B ₁	10006610
12	F6	Lauric Acid	10006626
12	F7	Palmitic Acid	10006627
12	F8	HEPC	10006695
12	F9	Heptanoyl thio-PC	10006809
12	F10	12-oxo Stearic Acid methyl ester	10007012
12	F11	CMPF	10007133
12	F12	<i>Unused</i>	
12	G1	<i>Unused</i>	
12	G2	N-Lignoceroyl Taurine	10007286
12	G3	N-Nervonoyl Taurine	10007287
12	G4	N-Docosanoyl Taurine	10007288
12	G5	<i>cis</i> -4,10,13,16-Docosatetraenoic Acid	10007289
12	G6	<i>cis</i> -7-Hexadecenoic Acid	10007290
12	G7	Arachidonoyl-N,N-dimethyl amide	10007293
12	G8	Arachidonoyl-N-methyl amide	10007294
12	G9	Arachidonoyl amide	10007295
12	G10	10-Thiastearic Acid	10007417
12	G11	Docosahexaenoyl Ethanolamide	10007534
12	G12	<i>Unused</i>	
12	H1	<i>Unused</i>	
12	H2	5 α -hydroxy-6-keto Cholesterol	10007601
12	H3	Δ^2 - <i>cis</i> Eicosenoic Acid	10007602
12	H4	N-Palmitoyl Dopamine	10007697
12	H5	N-(3-hydroxyphenyl)-Arachidonoyl amide	10007704
12	H6	PtdIns-(4)-P ₁ (1,2-dihexanoyl) (sodium salt)	10007757
12	H7	PtdIns-(3,4)-P ₂ (1,2-dihexanoyl) (sodium salt)	10007759
12	H8	PLC thio-PIP ₂ (sodium salt)	10007839
12	H9	8,11,14-Eicosatriynoic Acid	10007900
12	H10	Sphingosine (d20:1)	10007903
12	H11	Arachidonoyl-1-thio-Glycerol	10007904
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



Page 33 of 35

Plate	Well	Contents	Item Number
13	A1	<i>Unused</i>	
13	A2	17-hydroxy Heptadecanoic Acid	24650
13	A3	α -GalCer analog 8	26660
13	A4	Δ^7 -Avenasterol	30155
13	A5	Sphingosine (d18:1)	10007907
13	A6	20-carboxy Arachidonic Acid	10007912
13	A7	Sphinganine (d18:0)	10007945
13	A8	C16 Sphingomyelin (d18:1/16:0)	10007946
13	A9	D-erythro Lysosphingomyelin (d18:1)	10007947
13	A10	9-Nitrooleate	10008042
13	A11	10-Nitrooleate	10008043
13	A12	<i>Unused</i>	
13	B1	<i>Unused</i>	
13	B2	PtdIns-(5)-P ₁ (1,2-dihexanoyl) (sodium salt)	10008050
13	B3	Perifosine	10008112
13	B4	<i>cis</i> - Δ^2 -11-methyl-Dodecenoic Acid	10008123
13	B5	N-Cyclohexanecarbonyltetradecylamine	10008317
13	B6	<i>trans</i> - Δ^2 -11-methyl-Dodecenoic Acid	10008321
13	B7	all- <i>cis</i> -4,7,10,13,16-Docosapentaenoic Acid	10008335
13	B8	9(Z),11(E),13(E)-Octadecatrienoic Acid	10008349
13	B9	PtdIns-(3,4,5)-P ₃ (1,2-dihexanoyl) (ammonium salt)	10008390
13	B10	PtdIns-(3)-P ₁ (1,2-diocanoyl) (sodium salt)	10008394
13	B11	6 β -hydroxy Testosterone	10008519
13	B12	<i>Unused</i>	
13	C1	<i>Unused</i>	
13	C2	5- <i>cis</i> Iloprost	10008585
13	C3	Arachidonic Acid Leelamide	10008617
13	C4	Lauric Acid Leelamide	10008618
13	C5	oxy-Arachidonoyl Ethanolamide	10008642
13	C6	11 α -hydroxy Testosterone	10008647
13	C7	Pinolenic Acid	10008654
13	C8	13,14-dihydro-15-keto Prostaglandin J ₂	10008840
13	C9	1-Linoleoyl Glycerol	10008869
13	C10	N-Palmitoyl Glycine	10009020
13	C11	U-18666A	10009085
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 35

Plate	Well	Contents	Item Number
13	D1	<i>Unused</i>	
13	D2	CAY10509	10009167
13	D3	CAY10510	10009168
13	D4	(±)14(15)-EET-SI	10009286
13	D5	Ricinoleic Acid	10009369
13	D6	14(Z)-Eicosenoic Acid	10009375
13	D7	2-thio-Acetyl MAGE	10009651
13	D8	5(Z),8(Z),14(Z)-Eicosatrienoic Acid	10009733
13	D9	Palmitoleic Acid	10009871
13	D10	16-phenoxy Prostaglandin F _{2α} ethyl amide	10009875
13	D11	5(Z),11(Z),14(Z)-Eicosatrienoic Acid	10009999
13	D12	<i>Unused</i>	
13	E1	<i>Unused</i>	
13	E2	17-trifluoromethylphenyl trinor Prostaglandin F _{2α} ethyl amide	10010061
13	E3	17-trifluoromethylphenyl trinor Prostaglandin F _{2α} isopropyl ester	10010062
13	E4	11- <i>cis</i> Vaccenyl Acetate	10010101
13	E5	16-phenoxy tetranor Prostaglandin F _{2α} isopropyl ester	10010103
13	E6	15-deoxy-Δ ^{12,14} -Prostaglandin J ₂ -2-glyceryl ester	10010132
13	E7	1-Oleoyl-2-hydroxy- <i>sn</i> -glycero-3-PA	10010164
13	E8	12(Z)-Octadecenoic Acid methyl ester	10010180
13	E9	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PA	10010188
13	E10	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PA (sodium salt)	10010290
13	E11	Palmitoyl 3-carbacyclic Phosphatidic Acid	10010293
13	E12	<i>Unused</i>	
13	F1	<i>Unused</i>	
13	F2	Palmitoleoyl 3-carbacyclic Phosphatidic Acid	10010298
13	F3	Oleoyl 3-carbacyclic Phosphatidic Acid	10010299
13	F4	17-phenyl trinor Prostaglandin F _{2α} methyl amide	10010351
13	F5	Fluprostenol methyl amide	10010406
13	F6	13,14-dihydro-15-keto Prostaglandin D ₁	10010425
13	F7	AL 8810 ethyl amide	10010470
13	F8	(+)-Cloprostenol methyl amide	10010495
13	F9	Farnesyl Thiosalicylic Acid	10010501
13	F10	U-46619 Glycine methyl ester	10010522
13	F11	L- <i>threo</i> -Sphingosine (d18:1)	10010541
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



Page 35 of 35

Plate	Well	Contents	Item Number
13	G1	<i>Unused</i>	
13	G2	O-Arachidonoyl Glycidol	10010547
13	G3	16-phenoxy tetranor Prostaglandin F _{2α} methyl amide	10010562
13	G4	17-phenyl trinor Prostaglandin F _{2α} cyclopropyl amide	10010605
13	G5	17-phenoxy trinor Prostaglandin F _{2α} isopropyl ester	10010723
13	G6	17-phenoxy trinor Prostaglandin F _{2α} ethyl amide	10010742
13	G7	Capsaicin (technical grade)	10010743
13	G8	17-phenyl trinor Prostaglandin F _{2α} cyclopropyl methyl amide	10010810
13	G9	17-phenoxy trinor Prostaglandin F _{2α}	10010839
13	G10	Menaquinone 7	10011051
13	G11	Fulvestrant	10011269
13	G12	<i>Unused</i>	
13	H1	<i>Unused</i>	
13	H2	Chenodeoxycholic Acid	10011286
13	H3	Stearic Acid	10011298
13	H4	Glycerophospho-N-Arachidonoyl Ethanolamine	10011347
13	H5	Glycerophospho-N-Palmitoyl Ethanolamine	10011356
13	H6	Glycerophospho-N-Oleoyl Ethanolamine	10011357
13	H7	Phytomonic Acid	10012556
13	H8	7(Z),11(Z)-Heptacosadiene	10012567
13	H9	Glycerophospho-N-Eicosapentaenoyl Ethanolamine	90001842
13	H10	<i>Unused</i>	
13	H11	<i>Unused</i>	
13	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