

PRODUCT INFORMATION



Terpene Screening Library

Item No. 9003370 • Batch No. 0611615

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

Page 1 of 4

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Ursolic Acid	10072
1	A3	Forskolin	11018
1	A4	Betulin	11041
1	A5	Lupeol	11215
1	A6	Paxilline	11345
1	A7	β -acetyl-Boswellic Acid	11674
1	A8	Andrographolide	11679
1	A9	Bakuchiol	11684
1	A10	Betulinic Acid	11686
1	A11	β -Elemenic Acid	11712
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Oleanolic Acid	11726
1	B3	Neoandrographolide	11742
1	B4	Asiatic Acid	11818
1	B5	Madecassic Acid	11854
1	B6	Cafestol	13999
1	B7	Ingenol	14031
1	B8	Bilobalide	14272
1	B9	(-)-Huperzine A	14620
1	B10	Ginkgolide B	14636
1	B11	Cucurbitacin B	14820
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	Cucurbitacin E	14821
1	C3	Polygodial	14979
1	C4	Zerumbone	15400
1	C5	Ingenol-3-angelate	16207
1	C6	Ferutinin	16554
1	C7	Limonin	16932
1	C8	Phytol	17401
1	C9	Dehydrocostus lactone	18485
1	C10	β -Elemene	19641
1	C11	Juvenile Hormone III	19646
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, 04/21/2021

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 4

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Nomilin	20836
1	D3	β -Amyrin	20949
1	D4	Baccatin III	21404
1	D5	β -Caryophyllene	21572
1	D6	(-)-Caryophyllene oxide	21573
1	D7	α -Humulene	21574
1	D8	(\pm)-Linalool	21575
1	D9	α -Pinene	21576
1	D10	(-)- β -Pinene	21577
1	D11	Terpinolene	21712
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	(+)-Limonene	21719
1	E3	β -Myrcene	21720
1	E4	(-)-Perillyl Alcohol	21860
1	E5	Eucalyptol	22183
1	E6	Geranyl Acetate	22184
1	E7	Tubeimoside I	22276
1	E8	Incensole	22895
1	E9	<i>p</i> -Cymene	23159
1	E10	(+)-Isopulegol	23160
1	E11	(-)-Isopulegol	23161
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	(+)-Fenchol	23162
1	F3	(+)-3-Carene	23165
1	F4	(+)-Valencene	23167
1	F5	(-)-Guaiol	23172
1	F6	Terpineol	23173
1	F7	(1S)-(-)-Camphor	23176
1	F8	Isoborneol	23177
1	F9	α -Terpinene	23178
1	F10	α -Phellandrene	23179
1	F11	Cedrene	23180
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 4

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	Nerolidol	23182
1	G3	(+)-Menthol	23463
1	G4	(+)- β -Citronellol	23464
1	G5	(+)-Borneol	23467
1	G6	(-)-Borneol	23468
1	G7	γ -Terpinene	23722
1	G8	Columbin	24105
1	G9	(\pm)-Camphene	24478
1	G10	(+)-Nootkatone	24910
1	G11	Atractylenolide III	24911
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Abietic Acid	24927
1	H3	Ginkgolide A	24970
1	H4	Saikosaponin A	25026
1	H5	15,16-Dihydrotanshinone I	25051
1	H6	Atractylenolide I	25061
1	H7	Deacetylasperulosidic Acid	25100
1	H8	Oridonin	25665
1	H9	Benzoylaconine	25687
1	H10	Piperitenone	25752
1	H11	(\pm)-Menthol	25753
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	(+)-Cedrol	25769
2	A3	β -Eudesmol	25771
2	A4	(\pm)-Limonene	25773
2	A5	Nerol	25774
2	A6	Sabinene	25777
2	A7	Kirenol	26212
2	A8	Totarol	26412
2	A9	Curcumenol	26715
2	A10	Tanshinone I	26852
2	A11	Parthenolide	70080
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 4

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	Celastrol	70950
2	B3	Carnosic Acid	89820
2	B4	(-)- α -Bisabolol	9002495
2	B5	Leelamine	10006148
2	B6	Maslinic Acid	10009645
2	B7	<i>Unused</i>	
2	B8	<i>Unused</i>	
2	B9	<i>Unused</i>	
2	B10	<i>Unused</i>	
2	B11	<i>Unused</i>	
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	<i>Unused</i>	
2	C3	<i>Unused</i>	
2	C4	<i>Unused</i>	
2	C5	<i>Unused</i>	
2	C6	<i>Unused</i>	
2	C7	<i>Unused</i>	
2	C8	<i>Unused</i>	
2	C9	<i>Unused</i>	
2	C10	<i>Unused</i>	
2	C11	<i>Unused</i>	
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	<i>Unused</i>	
2	D3	<i>Unused</i>	
2	D4	<i>Unused</i>	
2	D5	<i>Unused</i>	
2	D6	<i>Unused</i>	
2	D7	<i>Unused</i>	
2	D8	<i>Unused</i>	
2	D9	<i>Unused</i>	
2	D10	<i>Unused</i>	
2	D11	<i>Unused</i>	
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