

PRODUCT INFORMATION



Fatty Acid Screening Library (96-Well)

Item No. 10504 • Batch No. 0574654

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

Page 1 of 3

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	N-(α -Linolenoyl) Tyrosine	10032
1	A3	N-Arachidonoyl-3-hydroxy- γ -Aminobutyric Acid	10158
1	A4	N-Oleoyl-L-Serine	13058
1	A5	Farnesyl Thiosalicylic Acid Amide	13175
1	A6	Myriocin	63150
1	A7	S-Farnesyl Thioacetic Acid	63260
1	A8	<i>cis</i> -Parinaric Acid	71430
1	A9	Traumatic Acid	88820
1	A10	Arachidonic Acid	90010
1	A11	Arachidonic Acid methyl ester	90014
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	N-Arachidonoyl-L-Alanine	90065
1	B3	N-Arachidonoyl- γ -Aminobutyric Acid	90067
1	B4	Eicosapentaenoic Acid	90110
1	B5	Eicosatetraenoic Acid	90120
1	B6	10(E),12(Z)-Conjugated Linoleic Acid	90145
1	B7	Linoleic Acid	90150
1	B8	Linoelaidic Acid	90160
1	B9	Docosapentaenoic Acid	90165
1	B10	Docosatrienoic Acid	90170
1	B11	13(Z)-Docosenoic Acid	90175
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	11(Z),14(Z),17(Z)-Eicosatrienoic Acid	90192
1	C3	α -Linolenic Acid	90210
1	C4	γ -Linolenic Acid	90220
1	C5	Dihomo- γ -Linolenic Acid	90230
1	C6	Dihomo- γ -Linolenic Acid ethyl ester	90236
1	C7	Elaidic Acid	90250
1	C8	Oleic Acid	90260
1	C9	17-Octadecynoic Acid	90270
1	C10	MEDICA 16	90290
1	C11	Adrenic Acid	90300
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, 04/27/2021

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 3

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Docosahexaenoic Acid	90310
1	D3	Stearidonic Acid	90320
1	D4	11(Z),14(Z)-Eicosadienoic Acid	90330
1	D5	Phytanic Acid	90360
1	D6	9(E),11(E)-Conjugated Linoleic Acid	90370
1	D7	2-Fluoro Palmitic Acid	90380
1	D8	2-hydroxy Myristic Acid	90390
1	D9	8-methyl Nonanoic Acid	9000310
1	D10	Linoleoyl Glycine	9000326
1	D11	Docosahexaenoyl Glycine	9000328
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Arachidic Acid	9000339
1	E3	Docosahexaenoic Acid ethyl ester	9090310
1	E4	Stearidonic Acid methyl ester	10005000
1	E5	N-Arachidonoyl-L-Serine	10005455
1	E6	Dihomo- γ -Linolenic Acid methyl ester	10006580
1	E7	Arachidonic Acid (sodium salt)	10006607
1	E8	Lauric Acid	10006626
1	E9	Palmitic Acid	10006627
1	E10	Oleic Acid-2,6-diisopropylanilide	10006782
1	E11	Stearidonic Acid ethyl ester	10006856
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	<i>cis</i> -7-Hexadecenoic Acid methyl ester	10006864
1	F3	Docosahexaenoic Acid methyl ester	10006865
1	F4	<i>cis</i> -4,10,13,16-Docosatetraenoic Acid methyl ester	10006866
1	F5	<i>cis</i> -4,10,13,16-Docosatetraenoic Acid	10007289
1	F6	<i>cis</i> -7-Hexadecenoic Acid	10007290
1	F7	Palmitic Acid methyl ester	10007358
1	F8	10-Thiastearic Acid	1007417
1	F9	Δ^2 - <i>cis</i> Eicosenoic Acid	10007602
1	F10	Δ^2 - <i>trans</i> Eicosenoic Acid	10007622
1	F11	CUDA	10007923
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 3

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	AUDA	10007927
1	G3	Stearic Acid ethyl ester	10008196
1	G4	Myristic Acid ethyl ester	10008197
1	G5	Linoleic Acid ethyl ester	10008198
1	G6	Linolenic Acid ethyl ester	10008199
1	G7	Arachidonic Acid ethyl ester	10008200
1	G8	Oleic Acid ethyl ester	10008201
1	G9	Palmitic Acid ethyl ester	10008202
1	G10	Lauric Acid ethyl ester	10008203
1	G11	Palmitoleic Acid ethyl ester	10008204
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	9(Z),11(E),13(E)-Octadecatrienoic Acid methyl ester	10008333
1	H3	9(Z),11(E),13(E)-Octadecatrienoic Acid ethyl ester	10008350
1	H4	Myristoleic Acid methyl ester	10008581
1	H5	Pinolenic Acid	10008654
1	H6	Pinolenic Acid methyl ester	10008656
1	H7	Pinolenic Acid ethyl ester	10008658
1	H8	Ricinoleic Acid methyl ester	10009729
1	H9	Palmitoleic Acid	10009871
1	H10	<i>cis</i> -12-Octadecenoic Acid methyl ester	10010180
1	H11	Stearic Acid	10011298
1	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