

PRODUCT INFORMATION



LipidLaunch™ Ionizable and Cationic Lipid Screening Library (96-Well)

Item No. 43949 • Batch No. 0813187

Page 1 of 6

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

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	1,2-Dioleoyloxy-3-dimethylamino-propane	15109
1	A3	1,2-Dioleoyl-3-dimethylammonium-propane	25726
1	A4	SM-102	33474
1	A5	ALC-0315	34337
1	A6	DLin-KC2-DMA	34363
1	A7	DLin-MC3-DMA	34364
1	A8	93-O17O	34366
1	A9	93-O17S	34367
1	A10	Lipid 5	34372
1	A11	L-319	35051
1	A12	Unused	
1	B1	Unused	
1	B2	Lipid A6	35052
1	B3	Lipid 29	35337
1	B4	YSK05	35786
1	B5	306O ₁₁₀	36698
1	B6	C12-200	36699
1	B7	cKK-E12	36700
1	B8	DLin-DMA	36701
1	B9	ATX-100	36935
1	B10	NT1-O14B	37095
1	B11	Lipid C24	37122
1	B12	Unused	
1	C1	Unused	
1	C2	9A1P9	37276
1	C3	CL4H6	37279
1	C4	FTT5	37413
1	C5	DOG-IM4	37441
1	C6	80-O16B	37564
1	C7	OF-02	37652
1	C8	Lipid A9	37667
1	C9	L202	37841
1	C10	OF-Deg-Lin	37853
1	C11	OF-C4-Deg-Lin	37856
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, 11/17/2025

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 6

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	AA3-DLin	37903
1	D3	246C10	37907
1	D4	TT3	37909
1	D5	Lipid 2,2(8,8) 4C CH ₃	37910
1	D6	306-N16B	37927
1	D7	Lipid OA2 (hydrochloride)	37987
1	D8	AL-A12	38001
1	D9	1O14	38150
1	D10	cKK-E15	38211
1	D11	OC2-K3-E10	38243
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	YK-009	38282
1	E3	1,2-Dimyristoyl-3-dimethylammonium-propane	38310
1	E4	1,2-Dipalmitoyl-3-dimethylammonium-propane	38311
1	E5	CAP 2 (hydrochloride)	38315
1	E6	Lipid CL1	38320
1	E7	C13-112-tetra-tail	38329
1	E8	Lipid 222	38338
1	E9	Lipid A4	38351
1	E10	A12-Iso5-2DC18	38586
1	E11	ALC-0315 analogous-1	38591
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	503O ₁₃	38628
1	F3	PL1	38647
1	F4	AA-T3A-C12	38648
1	F5	Lipid Catechol	38665
1	F6	Lipid M	38775
1	F7	C12-SPM	38784
1	F8	RCB-4-8	38803
1	F9	(S)-C12-200	38885
1	F10	RM 133-3	38917
1	F11	RM 137-15	38918
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 6

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	C14-4	38942
1	G3	ATX-001	39037
1	G4	Lipid R6	39130
1	G5	Lipid III-45	39243
1	G6	244cis	39284
1	G7	C3-K2-E14	39323
1	G8	C12-113	39335
1	G9	11-A-M	39458
1	G10	Lipid C3	39516
1	G11	DLin-MC2-DMA	39731
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	DLin-MC4-DMA	39732
1	H3	Lipid 331	39771
1	H4	Lipid C2	39783
1	H5	ALC-0315 N-oxide	39797
1	H6	Lipid C-1	39975
1	H7	9322-O16B	40208
1	H8	C14-A1	40256
1	H9	E12CA1A3	40415
1	H10	IR-117-17	40420
1	H11	Lipid 88	40493
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	Lipid DIM1	40512
2	A3	4-O10b1	40603
2	A4	80-O18	40611
2	A5	11-10-8	40747
2	A6	SIL Lipid	40801
2	A7	C14-O2	40805
2	A8	Lipid 2-10	40811
2	A9	C9-200	40862
2	A10	C10-200	40863
2	A11	Lipid U-101	40902
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 6

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	Lipid 29d	40913
2	B3	IZ-Chol	40963
2	B4	H1L1A1B3	41199
2	B5	Lipid HTO12	41202
2	B6	L-369	41232
2	B7	Q ₁ -SM-102 (iodide)	41239
2	B8	Q ₁ -ALC-0315 (iodide)	41240
2	B9	Lipid 7-1	41304
2	B10	C12-TLRa	41338
2	B11	O12-D3-I3	41385
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	119-23	41386
2	C3	DMT7	41389
2	C4	DLin-MeOH	41432
2	C5	Lipid A	41708
2	C6	CP-LC-0729	41929
2	C7	C14-490	42004
2	C8	A28-C6B2	42017
2	C9	Si5-N14	42520
2	C10	THP1 Lipid	42585
2	C11	31hP	42706
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	514O _{6,10}	42917
2	D3	BCP-NC2-C12	42940
2	D4	Lipid TG4C	42941
2	D5	Lipid ND-O1	43010
2	D6	E10i-494	43084
2	D7	Lipid 20b	43121
2	D8	Lipid PPz-2R1	43160
2	D9	DMA4-H228	43807
2	D10	Lipid C16-C	44085
2	D11	CF3-2N6-UC18	44140
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 6

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	BP Lipid 229	44269
2	E3	Lipid TD5	44285
2	E4	BP Lipid 700	44388
2	E5	PPZ-A10	9004144
2	E6	1,2(R)-Dioleloxy-3-dimethylamino-propane	9004302
2	E7	Lipid 1 HG3	9004529
2	E8	<i>Unused</i>	
2	E9	<i>Unused</i>	
2	E10	<i>Unused</i>	
2	E11	<i>Unused</i>	
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	<i>Unused</i>	
2	F3	<i>Unused</i>	
2	F4	<i>Unused</i>	
2	F5	<i>Unused</i>	
2	F6	<i>Unused</i>	
2	F7	<i>Unused</i>	
2	F8	<i>Unused</i>	
2	F9	<i>Unused</i>	
2	F10	<i>Unused</i>	
2	F11	<i>Unused</i>	
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	<i>Unused</i>	
2	G3	<i>Unused</i>	
2	G4	<i>Unused</i>	
2	G5	<i>Unused</i>	
2	G6	<i>Unused</i>	
2	G7	<i>Unused</i>	
2	G8	<i>Unused</i>	
2	G9	<i>Unused</i>	
2	G10	<i>Unused</i>	
2	G11	<i>Unused</i>	
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 6

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	<i>Unused</i>	
2	H3	<i>Unused</i>	
2	H4	<i>Unused</i>	
2	H5	<i>Unused</i>	
2	H6	<i>Unused</i>	
2	H7	<i>Unused</i>	
2	H8	<i>Unused</i>	
2	H9	<i>Unused</i>	
2	H10	<i>Unused</i>	
2	H11	<i>Unused</i>	
2	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