

PRODUCT INFORMATION



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

Item No. 10507 • Batch No. 0593991

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

Page 1 of 8

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Prostaglandin A ₁ -biotin	10013
1	A3	MK-571	10029
1	A4	PGPC	10044
1	A5	6(S)-Lipoxin A ₄	10049
1	A6	(±)12-HpETE	10138
1	A7	15-deoxy-Δ ^{12,14} -Prostaglandin J ₂ -biotin	10141
1	A8	HPF	10159
1	A9	15(S)-HETE Ethanolamide	10169
1	A10	(±)14(15)-EpETE	10173
1	A11	(±)16(17)-EpDPA	10174
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	(±)19(20)-EpDPA	10175
1	B3	13- <i>epi</i> -12-oxo Phytodienoic Acid	10195
1	B4	19(R)-hydroxy Prostaglandin A ₂	10295
1	B5	Prostaglandin A ₃	10310
1	B6	Prostaglandin D ₃	12990
1	B7	2,3-dinor Prostaglandin E ₁	13120
1	B8	13,14-dihydro-19(R)-hydroxy Prostaglandin E ₁	13660
1	B9	11β-Misoprostol	13825
1	B10	19(R)-hydroxy Prostaglandin E ₁	13910
1	B11	15(R),19(R)-hydroxy Prostaglandin E ₁	13920
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	Prostaglandin E ₂ <i>p</i> -benzamido phenyl ester	14054
1	C3	8- <i>iso</i> -15-keto Prostaglandin E ₂	14390
1	C4	11β-Prostaglandin E ₂	14510
1	C5	16,16-dimethyl Prostaglandin E ₂ <i>p</i> -(<i>p</i> -acetamidobenzamido) phenyl ester	14753
1	C6	tetranor-PGEM	14840
1	C7	19(R)-hydroxy Prostaglandin E ₂	14910
1	C8	15(R),19(R)-hydroxy Prostaglandin E ₂	14920
1	C9	20-hydroxy Prostaglandin E ₂	14950
1	C10	Prostaglandin E ₃	14990
1	C11	2,3-dinor-6-keto Prostaglandin F _{1α} (sodium salt)	15120
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, 08/12/2021

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 8

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Δ^{17} -6-keto Prostaglandin F _{1α}	15230
1	D3	6,15-diketo-13,14-dihydro Prostaglandin F _{1α}	15270
1	D4	19(R)-hydroxy Prostaglandin F _{1α}	15910
1	D5	2,3-dinor-8-iso Prostaglandin F _{2α}	16290
1	D6	(\pm)5-iPF _{2α} -VI	16300
1	D7	8-iso-13,14-dihydro-15-keto Prostaglandin F _{2α}	16380
1	D8	8-iso-15-keto Prostaglandin F _{2α}	16390
1	D9	U-51605	16465
1	D10	2,3-dinor-11 β -Prostaglandin F _{2α}	16530
1	D11	11 β -13,14-dihydro-15-keto Prostaglandin F _{2α}	16540
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Fluprostenol isopropyl ester	16769
1	E3	5- <i>trans</i> Fluprostenol isopropyl ester	16777
1	E4	15(S)-Fluprostenol isopropyl ester	16788
1	E5	tetranor-PGFM	16840
1	E6	19(R)-hydroxy Prostaglandin F _{2α}	16910
1	E7	15(R),19(R)-hydroxy Prostaglandin F _{2α}	16920
1	E8	20-hydroxy Prostaglandin F _{2α}	16950
1	E9	Prostaglandin F _{3α}	16990
1	E10	8-iso Prostaglandin F _{3α}	16992
1	E11	17- <i>trans</i> Prostaglandin F _{3α}	16995
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	Prostaglandin H ₁	17015
1	F3	Prostaglandin H ₂	17020
1	F4	Carbaprostacyclin	18210
1	F5	13,14-dehydro-15-cyclohexyl Carbaprostacyclin	18212
1	F6	Iloprost	18215
1	F7	Ciprostene (calcium salt)	18216
1	F8	Prostaglandin I ₃ (sodium salt)	18300
1	F9	Δ^{12} -Prostaglandin J ₂	18550
1	F10	Prostaglandin K ₁	18800
1	F11	Prostaglandin K ₂	18900
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 8

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	Carbocyclic Thromboxane A ₂	19010
1	G3	I-SAP	19021
1	G4	11-dehydro Thromboxane B ₂	19500
1	G5	11-dehydro-2,3-dinor Thromboxane B ₂	19510
1	G6	I-BOP	19600
1	G7	Thromboxane B ₃	19990
1	G8	11-dehydro Thromboxane B ₃	19995
1	G9	Leukotriene A ₃ methyl ester	20009
1	G10	Leukotriene A ₄ methyl ester	20010
1	G11	Leukotriene B ₃	20109
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Leukotriene B ₄	20110
1	H3	Leukotriene B ₄ Ethanolamide	20112
1	H4	Leukotriene B ₄ dimethyl amide	20115
1	H5	12- <i>epi</i> Leukotriene B ₃	20134
1	H6	12- <i>epi</i> Leukotriene B ₄	20135
1	H7	20-carboxy Leukotriene B ₄	20180
1	H8	20-hydroxy Leukotriene B ₄	20190
1	H9	Leukotriene F ₄	20520
1	H10	(±)5-HEPE	32200
1	H11	<i>Unused</i>	
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	5(S)-HEPE	32210
2	A3	(±)8-HEPE	32340
2	A4	(±)9-HEPE	32400
2	A5	9(S)-HEPE	32410
2	A6	(±)11-HEPE	32500
2	A7	11(R)-HEPE	32505
2	A8	11(S)-HEPE	32510
2	A9	(±)12-HEPE	32540
2	A10	(±)15-HEPE	32700
2	A11	15(S)-HEPE	32710
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 8

Plate	Well	Contents	Item Number
2	B1	<i>Unused</i>	
2	B2	(±)7-HDHA	33300
2	B3	(±)8-HDHA	33350
2	B4	(±)10-HDHA	33400
2	B5	(±)11-HDHA	33450
2	B6	(±)13-HDHA	33500
2	B7	(±)17-HDHA	33650
2	B8	(±)20-HDHA	33750
2	B9	(±)5-HETE	34210
2	B10	(±)5-HETE lactone	34215
2	B11	(±)5-HETE methyl ester	34220
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	5(S)-HETE	34230
2	C3	5(S)-HETE lactone	34240
2	C4	5-OxoETE	34250
2	C5	(±)8-HETE	34340
2	C6	(±)9-HETE	34400
2	C7	9(R)-HETE	34405
2	C8	9(S)-HETE	34410
2	C9	(±)11-HETE	34500
2	C10	11(R)-HETE	34505
2	C11	11(S)-HETE	34510
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	(±)12-HETE	34550
2	D3	12(R)-HETE	34560
2	D4	tetranor-12(R)-HETE	34565
2	D5	<i>Unused</i>	
2	D6	(±)15-HETE	34700
2	D7	15(R)-HETE	34710
2	D8	15(S)-HETE	34720
2	D9	15-OxoETE	34730
2	D10	5(S),6(R)-DiHETE	35200
2	D11	5(S),6(S)-DiHETE	35210
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 8

Plate	Well	Contents	Item Number
2	E1	<i>Unused</i>	
2	E2	6- <i>trans</i> Leukotriene B ₄	35250
2	E3	5(S),15(S)-DiHETE	35280
2	E4	8(S),15(S)-DiHETE	35370
2	E5	8(S)-HETrE	36360
2	E6	15(S)-HETrE	36720
2	E7	(±)11-HEDE	37500
2	E8	11(R)-HEDE	37505
2	E9	11(S)-HEDE	37510
2	E10	(±)15-HEDE	37700
2	E11	15(R)-HEDE	37710
2	E12	<i>Unused</i>	
2	F1	<i>Unused</i>	
2	F2	15(S)-HEDE	37720
2	F3	15-OxoEDE	37730
2	F4	(±)9-HODE	38400
2	F5	9(R)-HODE	38405
2	F6	9(S)-HODE	38410
2	F7	9-OxoODE	38420
2	F8	(±)13-HODE	38600
2	F9	13(R)-HODE	38605
2	F10	13(S)-HODE	38610
2	F11	9(S)-HOTrE	39420
2	F12	<i>Unused</i>	
2	G1	<i>Unused</i>	
2	G2	13(S)-HOTrE(γ)	39610
2	G3	13(S)-HOTrE	39620
2	G4	15(S)-HpEPE	42710
2	G5	5(S)-HpETE	44230
2	G6	12(S)-HpETE	44570
2	G7	13(S)-HpOTrE(γ)	45210
2	G8	13(S)-HpOTrE	45220
2	G9	15(S)-HpEDE	47720
2	G10	Cholesteryl Linoleate Hydroperoxides	48001
2	G11	9(S)-HpODE	48410
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 8

Plate	Well	Contents	Item Number
2	H1	<i>Unused</i>	
2	H2	13(S)-HpODE	48610
2	H3	(±)5(6)-EET	50211
2	H4	(±)11(12)-EET	50511
2	H5	(±)14(15)-EET	50651
2	H6	(±)17(18)-EpETE	50861
2	H7	(±)5,6-DHET	51211
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>	
3	A1	<i>Unused</i>	
3	A2	(±)5(6)-DiHET lactone	51212
3	A3	(±)8(9)-DiHET	51351
3	A4	(±)11(12)-DiHET	51511
3	A5	(±)14(15)-DiHET	51651
3	A6	(±)9(10)-EpOME	52400
3	A7	(±)12(13)-EpOME	52450
3	A8	(±)9(10)-DiHOME	53400
3	A9	D- <i>myo</i> -Inositol-1,4,5-triphosphate (potassium salt)	60960
3	A10	D- <i>myo</i> -Inositol-1,3,4,5-tetraphosphate (sodium salt)	60980
3	A11	PtdIns-(4,5)-P ₂ (1,2-dioctanoyl) (sodium salt)	64910
3	A12	<i>Unused</i>	
3	B1	<i>Unused</i>	
3	B2	PtdIns-(3,4,5)-P ₃ (1,2-dipalmitoyl) (sodium salt)	64920
3	B3	PtdIns-(3)-P ₁ (1,2-dipalmitoyl) (ammonium salt)	64921
3	B4	PtdIns-(3,4)-P ₂ (1,2-dipalmitoyl) (sodium salt)	64922
3	B5	PtdIns-(4)-P ₁ (1,2-dipalmitoyl) (ammonium salt)	64923
3	B6	PtdIns-(4,5)-P ₂ (1,2-dipalmitoyl) (ammonium salt)	64924
3	B7	PtdIns-(5)-P ₁ (1,2-dipalmitoyl) (ammonium salt)	64925
3	B8	U-75302	70705
3	B9	Lactacystin	70980
3	B10	Calcitriol	71820
3	B11	Hyperforin	75650
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 8

Plate	Well	Contents	Item Number
3	C1	<i>Unused</i>	
3	C2	5,6-dehydro Arachidonic Acid	90020
3	C3	20-HETE	90030
3	C4	5(6)-EpETE methyl ester	90114
3	C5	(±)4(5)-EpDPA methyl ester	90314
3	C6	Lipoxin A ₄	90410
3	C7	11-dehydro Thromboxane B ₂ -d ₄	319500
3	C8	2,3-dinor Thromboxane B ₁	10006330
3	C9	Prostaglandin D ₂ -biotin	10006697
3	C10	Prostaglandin E ₂ -biotin	10006987
3	C11	Prostaglandin F _{2α} -biotin	10006988
3	C12	<i>Unused</i>	
3	D1	<i>Unused</i>	
3	D2	(±)14(15)-DiHETE	10006998
3	D3	(±)17(18)-DiHETE	10006999
3	D4	(±)19(20)-DiHDPA	10007001
3	D5	tetranor-12(S)-HETE	10007207
3	D6	20-carboxy Arachidonic Acid	10007912
3	D7	9-Nitrooleate	10008042
3	D8	10-Nitrooleate	10008043
3	D9	PtdIns-(3)-P ₁ (1,2-dioctanoyl) (sodium salt)	10008394
3	D10	<i>Unused</i>	
3	D11	D- <i>myo</i> -Inositol-2,4-diphosphate (sodium salt)	10008419
3	D12	<i>Unused</i>	
3	E1	<i>Unused</i>	
3	E2	D- <i>myo</i> -Inositol-3,4,5-triphosphate (sodium salt)	10008425
3	E3	D- <i>myo</i> -Inositol-4-phosphate (ammonium salt)	10008437
3	E4	D- <i>myo</i> -Inositol-1,4-diphosphate (sodium salt)	10008438
3	E5	D- <i>myo</i> -Inositol-1,5-diphosphate (sodium salt)	10008441
3	E6	D- <i>myo</i> -Inositol-1,3-diphosphate (sodium salt)	10008443
3	E7	D- <i>myo</i> -Inositol-1,2,3,4-tetraphosphate (sodium salt)	10008448
3	E8	D- <i>myo</i> -Inositol-1,2,3,5-tetraphosphate (sodium salt)	10008449
3	E9	D- <i>myo</i> -Inositol-1,2,3,6-tetraphosphate (sodium salt)	10008450
3	E10	<i>Unused</i>	
3	E11	D- <i>myo</i> -Inositol-1,2,4,5,6-pentaphosphate (sodium salt)	10008452
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 8

Plate	Well	Contents	Item Number
3	F1	<i>Unused</i>	
3	F2	D-myo-Inositol-2,3,4,5-tetraphosphate (ammonium salt)	10008456
3	F3	(±)5(6)-EET Ethanolamide	10008596
3	F4	(±)8(9)-EET Ethanolamide	10008597
3	F5	(±)11(12)-EET Ethanolamide	10008598
3	F6	(±)14(15)-EET Ethanolamide	10008599
3	F7	20-HETE Ethanolamide	10008602
3	F8	tetranor-PGDM-d ₆	10009039
3	F9	<i>Unused</i>	
3	F10	<i>Unused</i>	
3	F11	<i>Unused</i>	
3	F12	<i>Unused</i>	
3	G1	<i>Unused</i>	
3	G2	<i>Unused</i>	
3	G3	<i>Unused</i>	
3	G4	<i>Unused</i>	
3	G5	<i>Unused</i>	
3	G6	<i>Unused</i>	
3	G7	<i>Unused</i>	
3	G8	<i>Unused</i>	
3	G9	<i>Unused</i>	
3	G10	<i>Unused</i>	
3	G11	<i>Unused</i>	
3	G12	<i>Unused</i>	
3	H1	<i>Unused</i>	
3	H2	<i>Unused</i>	
3	H3	<i>Unused</i>	
3	H4	<i>Unused</i>	
3	H5	<i>Unused</i>	
3	H6	<i>Unused</i>	
3	H7	<i>Unused</i>	
3	H8	<i>Unused</i>	
3	H9	<i>Unused</i>	
3	H10	<i>Unused</i>	
3	H11	<i>Unused</i>	
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