

PRODUCT INFORMATION



Prostaglandin Screening Library I (96-Well)

Item No. 10501 • Batch No. 0502309

Page 1 of 3

The composition of the library will always vary somewhat depending upon our inventory.

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	Prostaglandin F _{2α} -1-glyceryl ester	10139
1	A3	8-iso Prostaglandin F _{1b}	15370
1	A4	Prostaglandin F _{2α} Ethanolamide	16013
1	A5	Prostaglandin F _{2α} 1,15-lactone	16120
1	A6	13,14-dihydro Prostaglandin F _{2α}	16660
1	A7	15(R)-15-methyl Prostaglandin F _{2α}	16730
1	A8	Fluprostenol	16768
1	A9	Latanoprost (free acid)	16811
1	A10	17-phenyl trinor Prostaglandin F _{2α} amide	16821
1	A11	20-ethyl Prostaglandin F _{2α}	16940
1	A12	Unused	
1	B1	Unused	
1	B2	Prostaglandin F _{2α} serinol amide	10194
1	B3	Prostaglandin F _{1b}	15410
1	B4	Prostaglandin F _{2α} Alcohol methyl ether	16014
1	B5	5-trans Prostaglandin F _{2α}	16210
1	B6	13,14-dihydro-15-keto Prostaglandin F _{2α}	16670
1	B7	15(R)-15-methyl Prostaglandin F _{2α} methyl ester	16734
1	B8	Fluprostenol isopropyl ester	16769
1	B9	Latanoprost	16812
1	B10	Latanoprost ethyl amide	16822
1	B11	Unused	
1	B12	Unused	
1	C1	Unused	
1	C2	Prostaglandin F _{1α}	15010
1	C3	13,14-dihydro-15-keto Prostaglandin F _{1α}	15670
1	C4	Prostaglandin F _{2α} ethyl amide	16016
1	C5	5-trans Prostaglandin F _{2α} (tromethamine salt)	16220
1	C6	13,14-dihydro-15-keto Prostaglandin F _{2α} isopropyl ester	16671
1	C7	15(R)-Prostaglandin F _{2α}	16740
1	C8	16-phenyl tetranor Prostaglandin F _{2α}	16770
1	C9	5-trans Latanoprost	16813
1	C10	17-phenyl trinor Prostaglandin F _{2α} diethyl amide	16823
1	C11	Unused	
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, 03/23/2017

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	Prostaglandin F _{1α} Alcohol	15020
1	D3	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{1α}	15895
1	D4	Prostaglandin F _{2α} (tromethamine salt)	16020
1	D5	8-iso Prostaglandin F _{2α}	16350
1	D6	Unoprostone	16680
1	D7	15(S)-15-methyl Prostaglandin F _{2α}	16743
1	D8	5- <i>trans</i> Fluprostenol isopropyl ester	16777
1	D9	15(R)-17-phenyl trinor Prostaglandin F _{2α}	16814
1	D10	17-phenyl trinor Prostaglandin F _{2α} isopropyl ester	16824
1	D11	<i>Unused</i>	
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	2,3-dinor-6-keto Prostaglandin F _{1α} (sodium salt)	15120
1	E3	15(R),19(R)-hydroxy Prostaglandin F _{1α}	15920
1	E4	Prostaglandin F _{2α} diethyl amide	16023
1	E5	8-iso Prostaglandin F _{2b}	16370
1	E6	Unoprostone isopropyl ester	16681
1	E7	15(S)-15-methyl Prostaglandin F _{2α} methyl ester	16744
1	E8	15-keto Fluprostenol isopropyl ester	16786
1	E9	15-keto Latanoprost (free acid)	16815
1	E10	15(R)-17-phenyl trinor Prostaglandin F _{2α} isopropyl ester	16825
1	E11	<i>Unused</i>	
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	D ¹⁷ -6-keto Prostaglandin F _{1α}	15230
1	F3	Prostaglandin F _{2α}	16010
1	F4	Prostaglandin F _{2α} isopropyl ester	16030
1	F5	Prostaglandin F _{2b}	16410
1	F6	Lumula	16685
1	F7	16,16-dimethyl Prostaglandin F _{2α}	16750
1	F8	15(S)-Fluprostenol	16787
1	F9	15-keto Latanoprost	16816
1	F10	tetranor-PGFM	16840
1	F11	<i>Unused</i>	
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	6,15-diketo-13,14-dihydro Prostaglandin F _{1α}	15270
1	G3	Prostaglandin F _{2α} methyl ester	16011
1	G4	Prostaglandin F _{2α} dimethyl amide	16032
1	G5	Prostaglandin F _{2b} (tromethamine salt)	16420
1	G6	15-cyclohexyl pentanor Prostaglandin F _{2α}	16710
1	G7	16-phenoxy tetranor Prostaglandin F _{2α}	16760
1	G8	15(S)-Fluprostenol isopropyl ester	16788
1	G9	15(S)-Latanoprost	16818
1	G10	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{2α}	16895
1	G11	<i>Unused</i>	
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	8-iso Prostaglandin F _{1α}	15350
1	H3	Prostaglandin F _{2α} Alcohol	16012
1	H4	Prostaglandin F _{2α} dimethyl amine	16033
1	H5	11b-Prostaglandin F _{2α} Ethanolamide	16522
1	H6	15-keto Prostaglandin F _{2α}	16720
1	H7	(+)-Cloprostenol	16765
1	H8	17-phenyl trinor Prostaglandin F _{2α}	16810
1	H9	17-phenyl trinor Prostaglandin F _{2α} ethyl amide	16820
1	H10	15(R),19(R)-hydroxy Prostaglandin F _{2α}	16920
1	H11	<i>Unused</i>	
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