

PRODUCT INFORMATION



Prostaglandin Screening Library I (96-Well)

Item No. 10501 • Batch No. 0609348

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

Page 1 of 6

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	Prostaglandin F _{2α} -1-glyceryl ester	10139
1	A3	Prostaglandin F _{2α} serinol amide	10194
1	A4	Tafluprost ethyl ester	11612
1	A5	13,14-dihydro-16,16-difluoro Prostaglandin F _{2α}	13609
1	A6	13,14-epoxy Fluprostenol isopropyl ester	13679
1	A7	Prostaglandin F _{1α}	15010
1	A8	Prostaglandin F _{1α} Alcohol	15020
1	A9	2,3-dinor-6-keto Prostaglandin F _{1α} (sodium salt)	15120
1	A10	3-methoxy Prostaglandin F _{1α}	15170
1	A11	6-keto Prostaglandin F _{1α}	15210
1	A12	Unused	
1	B1	Unused	
1	B2	6,15-diketo-13,14-dihydro Prostaglandin F _{1α}	15270
1	B3	8-iso Prostaglandin F _{1α}	15350
1	B4	8-iso Prostaglandin F _{1β}	15370
1	B5	Prostaglandin F _{1β}	15410
1	B6	11β-Prostaglandin F _{1β}	15450
1	B7	11-deoxy Prostaglandin F _{1β}	15510
1	B8	11-deoxy Prostaglandin F _{1α}	15512
1	B9	13,14-dihydro Prostaglandin F _{1α}	15610
1	B10	13,14-dihydro-15-keto Prostaglandin F _{1α}	15670
1	B11	15-keto Prostaglandin F _{1α}	15710
1	B12	Unused	
1	C1	Unused	
1	C2	17,20-dimethyl Prostaglandin F _{1α}	15790
1	C3	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{1α}	15895
1	C4	Prostaglandin F _{2α}	16010
1	C5	Prostaglandin F _{2α} methyl ester	16011
1	C6	Prostaglandin F _{2α} Alcohol	16012
1	C7	Prostaglandin F _{2α} Ethanolamide	16013
1	C8	Prostaglandin F _{2α} Alcohol methyl ether	16014
1	C9	Prostaglandin F _{2α} ethyl amide	16016
1	C10	Prostaglandin F _{2α} diethyl amide	16023
1	C11	Prostaglandin F _{2α} isopropyl ester	16030
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, 02/01/2024

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	Prostaglandin F _{2α} dimethyl amide	16032
1	D3	Prostaglandin F _{2α} dimethyl amine	16033
1	D4	5- <i>trans</i> Prostaglandin F _{2α}	16210
1	D5	8- <i>iso</i> Prostaglandin F _{2α}	16350
1	D6	8- <i>iso</i> Prostaglandin F _{2β}	16370
1	D7	8- <i>iso</i> -15(R)-Prostaglandin F _{2α}	16395
1	D8	Prostaglandin F _{2β}	16410
1	D9	16,16-dimethyl Prostaglandin F _{2β}	16470
1	D10	11-deoxy Prostaglandin F _{2α}	16500
1	D11	11-deoxy Prostaglandin F _{2β}	16510
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	11β-Prostaglandin F _{2α}	16520
1	E3	11β-Prostaglandin F _{2α} Ethanolamide	16522
1	E4	13,14-dihydro Prostaglandin F _{2α}	16660
1	E5	13,14-dihydro-15-keto Prostaglandin F _{2α}	16670
1	E6	15-keto Prostaglandin F _{2α}	16720
1	E7	15(R)-15-methyl Prostaglandin F _{2α}	16730
1	E8	15(R)-15-methyl Prostaglandin F _{2α} methyl ester	16734
1	E9	15(R)-Prostaglandin F _{2α}	16740
1	E10	15(S)-15-methyl Prostaglandin F _{2α}	16743
1	E11	15(S)-15-methyl Prostaglandin F _{2α} methyl ester	16744
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	16,16-dimethyl Prostaglandin F _{2α}	16750
1	F3	16-phenoxy tetranor Prostaglandin F _{2α}	16760
1	F4	(+)-Cloprostenol (sodium salt)	16766
1	F5	Fluprostenol	16768
1	F6	Fluprostenol isopropyl ester	16769
1	F7	5- <i>trans</i> Fluprostenol	16776
1	F8	5- <i>trans</i> Fluprostenol isopropyl ester	16777
1	F9	15-keto Fluprostenol	16785
1	F10	15-keto Fluprostenol isopropyl ester	16786
1	F11	15(S)-Fluprostenol	16787
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	15(S)-Fluprostenol isopropyl ester	16788
1	G3	17-phenyl trinor Prostaglandin F _{2α}	16810
1	G4	Latanoprost (free acid)	16811
1	G5	Latanoprost	16812
1	G6	5- <i>trans</i> Latanoprost	16813
1	G7	15(R)-17-phenyl trinor Prostaglandin F _{2α}	16814
1	G8	15-keto Latanoprost (free acid)	16815
1	G9	15-keto Latanoprost	16816
1	G10	15-keto-17-phenyl trinor Prostaglandin F _{2α}	16817
1	G11	15(S)-Latanoprost	16818
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	17-phenyl trinor Prostaglandin F _{2α} ethyl amide	16820
1	H3	17-phenyl trinor Prostaglandin F _{2α} amide	16821
1	H4	Latanoprost ethyl amide	16822
1	H5	17-phenyl trinor Prostaglandin F _{2α} diethyl amide	16823
1	H6	17-phenyl trinor Prostaglandin F _{2α} isopropyl ester	16824
1	H7	15(R)-17-phenyl trinor Prostaglandin F _{2α} isopropyl ester	16825
1	H8	17-trifluoromethylphenyl trinor Prostaglandin F _{2α}	16890
1	H9	17-trifluoromethylphenyl-13,14-dihydro trinor Prostaglandin F _{2α}	16895
1	H10	20-ethyl Prostaglandin F _{2α}	16940
1	H11	15-keto Cloprostenol isopropyl ester	24929
1	H12	<i>Unused</i>	
2	A1	<i>Unused</i>	
2	A2	(+)-5- <i>trans</i> Cloprostenol isopropyl ester	24930
2	A3	17-phenyl trinor Prostaglandin F _{2α} dimethyl amide	25241
2	A4	17-phenyl trinor Prostaglandin F _{2α} isopropyl amide	25242
2	A5	Latanoprost amide	25245
2	A6	Latanoprost dimethyl amide	25247
2	A7	1a-monohomo-17-phenyl trinor Prostaglandin F _{2α}	25392
2	A8	17-phenyl trinor Prostaglandin F _{2α} glycnamide	25470
2	A9	17-phenyl trinor Prostaglandin F _{2α} glycnamide methyl ester	25471
2	A10	15(R)-Prostaglandin F _{1α}	29549
2	A11	17-phenyl trinor Prostaglandin F _{2α} cyclohexyl amide	9000686
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	Tafluprost ethyl amide	9000843
2	B3	Fluprostenol serinol amide	10004236
2	B4	17-phenyl trinor Prostaglandin F _{2α} serinol amide	10004237
2	B5	(+)-5- <i>trans</i> Cloprostenol	10004970
2	B6	Tafluprost (free acid)	10005439
2	B7	Tafluprost	10005440
2	B8	8- <i>iso</i> Prostaglandin F _{2α} Ethanolamide	10005764
2	B9	<i>ent</i> -Prostaglandin F _{2α}	10008122
2	B10	15(R)-17-phenyl trinor Prostaglandin F _{2α} ethyl amide	10008127
2	B11	5- <i>trans</i> -17-phenyl trinor Prostaglandin F _{2α} ethyl amide	10008132
2	B12	<i>Unused</i>	
2	C1	<i>Unused</i>	
2	C2	8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2α}	10008435
2	C3	8- <i>iso</i> -17-phenyl trinor Prostaglandin F _{2β}	10008436
2	C4	8- <i>iso</i> -15-keto Prostaglandin F _{2β}	10008539
2	C5	16-phenoxy Prostaglandin F _{2α} ethyl amide	10009875
2	C6	(+)-Cloprostenol isopropyl ester	10010016
2	C7	17-trifluoromethylphenyl trinor Prostaglandin F _{2α} ethyl amide	10010061
2	C8	17-trifluoromethylphenyl trinor Prostaglandin F _{2α} isopropyl ester	10010062
2	C9	16-phenoxy tetranor Prostaglandin F _{2α} methyl ester	10010102
2	C10	16-phenoxy tetranor Prostaglandin F _{2α} isopropyl ester	10010103
2	C11	17-phenyl trinor Prostaglandin F _{2α} methyl ester	10010110
2	C12	<i>Unused</i>	
2	D1	<i>Unused</i>	
2	D2	17-trifluoromethylphenyl trinor Prostaglandin F _{2α} methyl ester	10010111
2	D3	(+)-Cloprostenol methyl ester	10010115
2	D4	Fluprostenol methyl ester	10010151
2	D5	17-phenyl trinor Prostaglandin F _{2α} methyl amide	10010351
2	D6	15-keto-17-phenyl trinor Prostaglandin F _{2α} ethyl amide	10010405
2	D7	Fluprostenol methyl amide	10010406
2	D8	(+)-Cloprostenol methyl amide	10010495
2	D9	16-phenoxy tetranor Prostaglandin F _{2α} methyl amide	10010562
2	D10	17-phenyl trinor Prostaglandin F _{2α} cyclopropyl amide	10010605
2	D11	17-phenoxy trinor Prostaglandin F _{2α} isopropyl ester	10010723
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	17-phenoxy trinor Prostaglandin F _{2α} ethyl amide	10010742
2	E3	17-phenyl trinor Prostaglandin F _{2α} cyclopropyl methyl amide	10010810
2	E4	17-phenoxy trinor Prostaglandin F _{2α}	10010839
2	E5	<i>Unused</i>	
2	E6	<i>Unused</i>	
2	E7	<i>Unused</i>	
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	Unused	
2	H2	Unused	
2	H3	Unused	
2	H4	Unused	
2	H5	Unused	
2	H6	Unused	
2	H7	Unused	
2	H8	Unused	
2	H9	Unused	
2	H10	Unused	
2	H11	Unused	
2	H12	Unused	

*Primary Target is based on available IC_{50} and/or K_i values cited in the literature

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