

Product Information



Prostaglandin Screening Library II (96-Well)

Item No. 10502 • Batch No. 0469450

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

Page 1 of 3

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	15(R)-Prostaglandin D ₂	10118
1	A3	Prostaglandin D ₂ -1-glyceryl ester	12015
1	A4	Prostaglandin E ₁ Ethanolamide	13012
1	A5	15-keto Prostaglandin E ₁	13680
1	A6	Prostaglandin E ₂	14010
1	A7	11β-Prostaglandin E ₂	14510
1	A8	Sulprostone	14765
1	A9	11-keto Fluprostenol	16783
1	A10	<i>Unused</i>	
1	A11	<i>Unused</i>	
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	Prostaglandin E ₂ -1-glyceryl ester	10140
1	B3	11-deoxy-11-methylene Prostaglandin D ₂	12410
1	B4	Prostaglandin E ₁ Alcohol	13020
1	B5	15(S)-15-methyl Prostaglandin E ₁	13730
1	B6	Prostaglandin E ₂ methyl ester	14011
1	B7	13,14-dihydro-15-keto Prostaglandin E ₂	14650
1	B8	16-phenyl tetranor Prostaglandin E ₂	14770
1	B9	(R)-Butaprost (free acid)	10006045
1	B10	<i>Unused</i>	
1	B11	<i>Unused</i>	
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	Prostaglandin D ₂ serinol amide	10192
1	C3	13,14-dihydro-15-keto Prostaglandin D ₂	12610
1	C4	1a,1b-dihomo Prostaglandin E ₁	13050
1	C5	(R)-Butaprost	13745
1	C6	Prostaglandin E ₂ Ethanolamide	14012
1	C7	15(R)-Prostaglandin E ₂	14710
1	C8	17-phenyl trinor Prostaglandin E ₂	14810
1	C9	Prostaglandin E ₂ -biotin	10006987
1	C10	<i>Unused</i>	
1	C11	<i>Unused</i>	
1	C12	<i>Unused</i>	

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/16/2015

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

Product Information

Page 2 of 3



Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Prostaglandin E ₂ serinol amide	10193
1	D3	Δ^{12} -Prostaglandin D ₂	12650
1	D4	6-keto Prostaglandin E ₁	13260
1	D5	CAY10408	13747
1	D6	Prostaglandin E ₂ <i>p</i> -acetamidophenyl ester	14053
1	D7	15-keto Prostaglandin E ₂	14720
1	D8	tetranor-PGEM	14840
1	D9	<i>ent</i> -Prostaglandin E ₂	10008294
1	D10	<i>Unused</i>	
1	D11	<i>Unused</i>	
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	Prostaglandin D ₁	12000
1	E3	15(R)-15-methyl Prostaglandin D ₂	12720
1	E4	8- <i>iso</i> Prostaglandin E ₁	13360
1	E5	16,16-dimethyl Prostaglandin E ₁	13750
1	E6	Prostaglandin E ₂ <i>p</i> -benzamido-phenyl ester	14054
1	E7	15(R)-15-methyl Prostaglandin E ₂	14725
1	E8	19(R)-hydroxy Prostaglandin E ₂	14910
1	E9	Prostaglandin D ₂ methyl ester	10008385
1	E10	<i>Unused</i>	
1	E11	<i>Unused</i>	
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	Prostaglandin D ₁ Alcohol	12002
1	F3	15(S)-15-methyl Prostaglandin D ₂	12730
1	F4	13,14-dihydro Prostaglandin E ₁	13610
1	F5	16-phenyl tetranor Prostaglandin E ₁	13770
1	F6	5- <i>trans</i> Prostaglandin E ₂	14210
1	F7	15(S)-15-methyl Prostaglandin E ₂	14730
1	F8	20-ethyl Prostaglandin E ₂	14940
1	F9	3-methoxy Limaprost	10008445
1	F10	<i>Unused</i>	
1	F11	<i>Unused</i>	
1	F12	<i>Unused</i>	

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

Product Information

Page 3 of 3



Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	Prostaglandin D ₂	12010
1	G3	17-phenyl trinor Prostaglandin D ₂	12810
1	G4	13,14-dihydro-15(R)-Prostaglandin E ₁	13630
1	G5	Limaprost	13810
1	G6	8- <i>iso</i> Prostaglandin E ₂	14350
1	G7	16,16-dimethyl Prostaglandin E ₂	14750
1	G8	Prostaglandin E ₃	14990
1	G9	8- <i>iso</i> -16-cyclohexyl-tetranor Prostaglandin E ₂	10009278
1	G10	<i>Unused</i>	
1	G11	<i>Unused</i>	
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	Prostaglandin D ₂ Ethanolamide	12012
1	H3	Prostaglandin E ₁	13010
1	H4	13,14-dihydro-15-keto Prostaglandin E ₁	13650
1	H5	Misoprostol	13820
1	H6	8- <i>iso</i> Prostaglandin E ₂ isopropyl ester	14352
1	H7	16,16-dimethyl Prostaglandin E ₂ <i>p</i> -(<i>p</i> -acetamidobenzamido) phenyl ester	14753
1	H8	17- <i>trans</i> Prostaglandin E ₃	14995
1	H9	13,14-dihydro-15-keto Prostaglandin D ₁	10010425
1	H10	<i>Unused</i>	
1	H11	<i>Unused</i>	
1	H12	<i>Unused</i>	

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com