

# PRODUCT INFORMATION



## Prostaglandin Screening Library II (96-Well)

Item No. 10502 • Batch No. 0584638

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

Page 1 of 2

Plate	Well	Contents	Item Number
1	A1	Unused	
1	A2	Prostaglandin E <sub>2</sub> -1-glycerol ester	10140
1	A3	Prostaglandin E <sub>2</sub> serinol amide	10193
1	A4	Prostaglandin D <sub>1</sub>	12000
1	A5	Prostaglandin D <sub>1</sub> Alcohol	12002
1	A6	Prostaglandin D <sub>2</sub>	12010
1	A7	Prostaglandin D <sub>2</sub> Ethanolamide	12012
1	A8	Prostaglandin D <sub>2</sub> -1-glycerol ester	12015
1	A9	11-deoxy-11-methylene Prostaglandin D <sub>2</sub>	12410
1	A10	13,14-dihydro-15-keto Prostaglandin D <sub>2</sub>	12610
1	A11	Δ <sup>12</sup> -Prostaglandin D <sub>2</sub>	12650
1	A12	Unused	
1	B1	Unused	
1	B2	15(S)-15-methyl Prostaglandin D <sub>2</sub>	12730
1	B3	17-phenyl trinor Prostaglandin D <sub>2</sub>	12810
1	B4	Prostaglandin E <sub>1</sub>	13010
1	B5	Prostaglandin E <sub>1</sub> Ethanolamide	13012
1	B6	Prostaglandin E <sub>1</sub> Alcohol	13020
1	B7	1a,1b-dihomo Prostaglandin E <sub>1</sub>	13050
1	B8	8-iso Prostaglandin E <sub>1</sub>	13360
1	B9	13,14-dihydro Prostaglandin E <sub>1</sub>	13610
1	B10	13,14-dihydro-15(R)-Prostaglandin E <sub>1</sub>	13630
1	B11	13,14-dihydro-15-keto Prostaglandin E <sub>1</sub>	13650
1	B12	Unused	
1	C1	Unused	
1	C2	15-keto Prostaglandin E <sub>1</sub>	13680
1	C3	15(S)-15-methyl Prostaglandin E <sub>1</sub>	13730
1	C4	(R)-Butaprost	13745
1	C5	CAY10408	13747
1	C6	16,16-dimethyl Prostaglandin E <sub>1</sub>	13750
1	C7	16-phenyl tetranor Prostaglandin E <sub>1</sub>	13770
1	C8	Limaprost	13810
1	C9	Misoprostol	13820
1	C10	Prostaglandin E <sub>2</sub>	14010
1	C11	Prostaglandin E <sub>2</sub> methyl ester	14011
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, 04/21/2021

#### 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 2

Plate	Well	Contents	Item Number
1	D1	<i>Unused</i>	
1	D2	Prostaglandin E <sub>2</sub> Ethanolamide	14012
1	D3	Prostaglandin E <sub>2</sub> <i>p</i> -acetamidophenyl ester	14053
1	D4	Prostaglandin E <sub>2</sub> <i>p</i> -benzamidophenyl ester	14054
1	D5	5- <i>trans</i> Prostaglandin E <sub>2</sub>	14210
1	D6	8- <i>iso</i> Prostaglandin E <sub>2</sub>	14350
1	D7	8- <i>iso</i> Prostaglandin E <sub>2</sub> isopropyl ester	14352
1	D8	11 $\beta$ -Prostaglandin E <sub>2</sub>	14510
1	D9	13,14-dihydro-15-keto Prostaglandin E <sub>2</sub>	14650
1	D10	15(R)-Prostaglandin E <sub>2</sub>	14710
1	D11	15-keto Prostaglandin E <sub>2</sub>	14720
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	15(R)-15-methyl Prostaglandin E <sub>2</sub>	14725
1	E3	15(S)-15-methyl Prostaglandin E <sub>2</sub>	14730
1	E4	16,16-dimethyl Prostaglandin E <sub>2</sub>	14750
1	E5	Sulprostone	14765
1	E6	17-phenyl trinor Prostaglandin E <sub>2</sub>	14810
1	E7	tetranor-PGEM	14840
1	E8	19(R)-hydroxy Prostaglandin E <sub>2</sub>	14910
1	E9	20-ethyl Prostaglandin E <sub>2</sub>	14940
1	E10	Prostaglandin E <sub>3</sub>	14990
1	E11	17- <i>trans</i> Prostaglandin E <sub>3</sub>	14995
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	11-keto Fluprostenol	16783
1	F3	(R)-Butaprost (free acid)	10006045
1	F4	Prostaglandin E <sub>2</sub> -biotin	10006987
1	F5	ent-Prostaglandin E <sub>2</sub>	10008294
1	F6	3-methoxy Limaprost	10008445
1	F7	8- <i>iso</i> -16-cyclohexyl-tetranor Prostaglandin E <sub>2</sub>	10009278
1	F8	13,14-dihydro-15-keto Prostaglandin D <sub>1</sub>	10010425
1	F9	<i>Unused</i>	
1	F10	<i>Unused</i>	
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