

# PRODUCT INFORMATION



## Prostaglandin Screening Library III (96-Well)

Item No. 10503 • Batch No. 0625990

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

Page 1 of 3

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	Prostaglandin A <sub>1</sub>	10010
1	A3	8-iso Prostaglandin A <sub>1</sub>	10035
1	A4	15-deoxy- $\Delta^{12,14}$ -Prostaglandin A <sub>1</sub>	10065
1	A5	15- <i>epi</i> Prostaglandin A <sub>1</sub>	10070
1	A6	16,16-dimethyl Prostaglandin A <sub>1</sub>	10080
1	A7	Prostaglandin A <sub>2</sub>	10210
1	A8	8-iso Prostaglandin A <sub>2</sub>	10235
1	A9	13,14-dihydro-15-keto Prostaglandin A <sub>2</sub>	10260
1	A10	15-deoxy- $\Delta^{12,14}$ -Prostaglandin A <sub>2</sub>	10265
1	A11	15(R)-15-methyl Prostaglandin A <sub>2</sub>	10270
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	16,16-dimethyl Prostaglandin A <sub>2</sub>	10280
1	B3	16-phenoxy tetranor Prostaglandin A <sub>2</sub>	10285
1	B4	17-phenyl trinor Prostaglandin A <sub>2</sub>	10288
1	B5	17-phenyl trinor-13,14-dihydro Prostaglandin A <sub>2</sub>	10290
1	B6	19(R)-hydroxy Prostaglandin A <sub>2</sub>	10295
1	B7	Prostaglandin B <sub>1</sub>	11110
1	B8	Prostaglandin B <sub>2</sub>	11210
1	B9	19(R)-hydroxy Prostaglandin B <sub>2</sub>	11910
1	B10	Treprostinil (diethanolamine salt)	11927
1	B11	BW 245C	12050
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	BW 246C	12055
1	C3	Bicyclo Prostaglandin E <sub>1</sub>	13530
1	C4	17-phenyl trinor Prostaglandin F <sub>2<math>\alpha</math></sub> 1,15-lactone	13992
1	C5	Bicyclo Prostaglandin E <sub>2</sub>	14530
1	C6	Prostaglandin F <sub>2<math>\alpha</math></sub> 1,11-lactone	16110
1	C7	Prostaglandin F <sub>2<math>\alpha</math></sub> 1,15-lactone	16120
1	C8	Prostaglandin F <sub>2<math>\alpha</math></sub> 1,9-lactone	16130
1	C9	5- <i>trans</i> U-44069	16442
1	C10	5- <i>trans</i> U-46619	16452
1	C11	6 $\alpha$ -Prostaglandin I <sub>1</sub>	18110
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, 06/19/2025

#### 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	6 $\beta$ -Prostaglandin I <sub>1</sub>	18120
1	D3	Carbaprostacyclin	18210
1	D4	5- <i>cis</i> Carbaprostacyclin	18211
1	D5	Iloprost	18215
1	D6	Ciprostene (calcium salt)	18216
1	D7	Prostaglandin I <sub>2</sub> (sodium salt)	18220
1	D8	Piriprost (potassium salt)	18228
1	D9	Beraprost (sodium salt)	18230
1	D10	Prostaglandin J <sub>2</sub>	18500
1	D11	$\Delta^{12}$ -Prostaglandin J <sub>2</sub>	18550
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	15-deoxy- $\Delta^{12,14}$ -Prostaglandin J <sub>2</sub>	18570
1	E3	15-deoxy- $\Delta^{12,14}$ -Prostaglandin J <sub>2</sub>	18570.1
1	E4	15-deoxy- $\Delta^{12,14}$ -Prostaglandin J <sub>2</sub> Glutathione	18580
1	E5	CAY10410	18590
1	E6	Pinane Thromboxane A <sub>2</sub>	19020
1	E7	SQ 29,548	19025
1	E8	I-SAP	19021
1	E9	Thromboxane B <sub>2</sub>	19030
1	E10	11-dehydro Thromboxane B <sub>2</sub>	19500
1	E11	15(R)-Prostaglandin I <sub>2</sub> (sodium salt)	21586
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	15(S)-acetate Prostaglandin A <sub>2</sub> methyl ester	28195
1	F3	15-keto Prostaglandin A <sub>1</sub>	9000142
1	F4	A <sub>1</sub> -Phytoprostane-I	9000593
1	F5	Thromboxane B <sub>1</sub>	10006610
1	F6	15(R)-Pinane Thromboxane A <sub>2</sub>	10008510
1	F7	5- <i>cis</i> Iloprost	10008585
1	F8	5- <i>cis</i> -15(R)-Iloprost	10008588
1	F9	15-keto Iloprost	10008616
1	F10	13,14-dihydro-15-keto Prostaglandin J <sub>2</sub>	10008840
1	F11	Taprostene (free acid)	10011348
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	Unused	
1	G2	Unused	
1	G3	Unused	
1	G4	Unused	
1	G5	Unused	
1	G6	Unused	
1	G7	Unused	
1	G8	Unused	
1	G9	Unused	
1	G10	Unused	
1	G11	Unused	
1	G12	Unused	
1	H1	Unused	
1	H2	Unused	
1	H3	Unused	
1	H4	Unused	
1	H5	Unused	
1	H6	Unused	
1	H7	Unused	
1	H8	Unused	
1	H9	Unused	
1	H10	Unused	
1	H11	Unused	
1	H12	Unused	

## 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