

PRODUCT INFORMATION



D-myo-Inositol Phosphate Screening Library (96-Well)

Item No. 9003895 • Batch No. 0624645

This library contains 5 nmol of lyophilized powder, sufficient for 0.1 mM stock solutions in 50 µl of an aqueous solution of choice.

Page 1 of 3

Plate	Well	Contents	Item Number
1	A1	<i>Unused</i>	
1	A2	D-myo-Inositol-1,4,5-triphosphate (potassium salt)	60960
1	A3	D-myo-Inositol-1,3,4-triphosphate (sodium salt)	60972
1	A4	D-myo-Inositol-1,3,4,5-tetraphosphate (sodium salt)	60980
1	A5	D-myo-Inositol-1-phosphate (cyclohexyl ammonium salt)	9001248
1	A6	D-myo-Inositol-1-phosphate (sodium salt)	10007777
1	A7	D-myo-Inositol-3-phosphate (sodium salt)	10007778
1	A8	D-myo-Inositol-2,4,5-triphosphate (sodium salt)	10007779
1	A9	D-myo-Inositol-1,2,6-triphosphate (sodium salt)	10007780
1	A10	D-myo-Inositol-1,3,5-triphosphate (sodium salt)	10007781
1	A11	D-myo-Inositol-3,4,5,6-tetraphosphate (sodium salt)	10007782
1	A12	<i>Unused</i>	
1	B1	<i>Unused</i>	
1	B2	D-myo-Inositol-1,4,5,6-tetraphosphate (sodium salt)	10007783
1	B3	D-myo-Inositol-1,3,4,5,6-pentaphosphate (sodium salt)	10007784
1	B4	D-myo-Inositol-1,4,5-triphosphate (sodium salt)	10008205
1	B5	D-myo-Inositol-1,2,3,4,5,6-hexaphosphate (sodium salt)	10008415
1	B6	D-myo-Inositol-4,5-diphosphate (sodium salt)	10008418
1	B7	D-myo-Inositol-2,4-diphosphate (sodium salt)	10008419
1	B8	D-myo-Inositol-1,5,6-triphosphate (sodium salt)	10008422
1	B9	D-myo-Inositol-2,3,5-triphosphate (ammonium salt)	10008423
1	B10	D-myo-Inositol-2,5,6-triphosphate (sodium salt)	10008424
1	B11	D-myo-Inositol-3,4,5-triphosphate (sodium salt)	10008425
1	B12	<i>Unused</i>	
1	C1	<i>Unused</i>	
1	C2	D-myo-Inositol-1,4,6-triphosphate (sodium salt)	10008427
1	C3	D-myo-Inositol-4-phosphate (ammonium salt)	10008437
1	C4	D-myo-Inositol-1,4-diphosphate (sodium salt)	10008438
1	C5	D-myo-Inositol-1,2-diphosphate (sodium salt)	10008439
1	C6	D-myo-Inositol-1,5-diphosphate (sodium salt)	10008441
1	C7	D-myo-Inositol-1,3,4,6-tetraphosphate (ammonium salt)	10008442
1	C8	D-myo-Inositol-1,3-diphosphate (sodium salt)	10008443
1	C9	D-myo-Inositol-1,2,5,6-tetraphosphate (sodium salt)	10008444
1	C10	D-myo-Inositol-1,2,3,4-tetraphosphate (sodium salt)	10008448
1	C11	D-myo-Inositol-1,2,3,5-tetraphosphate (sodium salt)	10008449
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, 04/11/2025

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	D-myo-Inositol-1,2,3,6-tetraphosphate (sodium salt)	10008450
1	D3	D-myo-Inositol-1,2,4,5-tetraphosphate (sodium salt)	10008451
1	D4	D-myo-Inositol-1,2,4,5,6-pentaphosphate (sodium salt)	10008452
1	D5	D-myo-Inositol-1,2,3,5,6-pentaphosphate (sodium salt)	10008453
1	D6	D-myo-Inositol-2,3,5,6-tetraphosphate (sodium salt)	10008455
1	D7	D-myo-Inositol-2,3,4,5-tetraphosphate (ammonium salt)	10008456
1	D8	D-myo-Inositol-1,3,4,5,6-pentaphosphate (ammonium salt)	10009851
1	D9	<i>Unused</i>	
1	D10	<i>Unused</i>	
1	D11	<i>Unused</i>	
1	D12	<i>Unused</i>	
1	E1	<i>Unused</i>	
1	E2	<i>Unused</i>	
1	E3	<i>Unused</i>	
1	E4	<i>Unused</i>	
1	E5	<i>Unused</i>	
1	E6	<i>Unused</i>	
1	E7	<i>Unused</i>	
1	E8	<i>Unused</i>	
1	E9	<i>Unused</i>	
1	E10	<i>Unused</i>	
1	E11	<i>Unused</i>	
1	E12	<i>Unused</i>	
1	F1	<i>Unused</i>	
1	F2	<i>Unused</i>	
1	F3	<i>Unused</i>	
1	F4	<i>Unused</i>	
1	F5	<i>Unused</i>	
1	F6	<i>Unused</i>	
1	F7	<i>Unused</i>	
1	F8	<i>Unused</i>	
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

PRODUCT INFORMATION



Page 3 of 3

Plate	Well	Contents	Item Number
1	G1	<i>Unused</i>	
1	G2	<i>Unused</i>	
1	G3	<i>Unused</i>	
1	G4	<i>Unused</i>	
1	G5	<i>Unused</i>	
1	G6	<i>Unused</i>	
1	G7	<i>Unused</i>	
1	G8	<i>Unused</i>	
1	G9	<i>Unused</i>	
1	G10	<i>Unused</i>	
1	G11	<i>Unused</i>	
1	G12	<i>Unused</i>	
1	H1	<i>Unused</i>	
1	H2	<i>Unused</i>	
1	H3	<i>Unused</i>	
1	H4	<i>Unused</i>	
1	H5	<i>Unused</i>	
1	H6	<i>Unused</i>	
1	H7	<i>Unused</i>	
1	H8	<i>Unused</i>	
1	H9	<i>Unused</i>	
1	H10	<i>Unused</i>	
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