

D-myo-Inositol 1-monophosphate . dipotassium salt

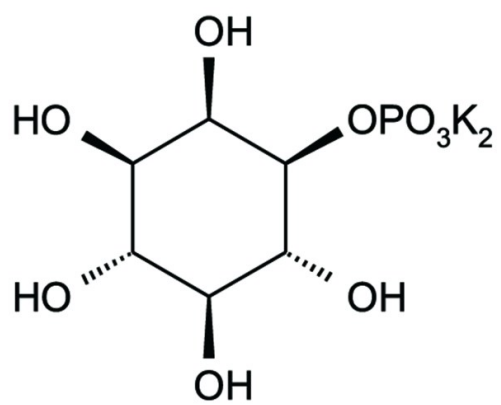
Ordering Information

[Order Online »](#)

ALX-307-026-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and moisture. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ins(1)P1 . 2K
Appearance	White to off-white solid.
CAS	313263-16-0 / 15421-51-9 (H)
Formula	$C_6H_{11}O_9P \cdot 2K$
Identity	Determined by 1H -NMR, ^{31}P -NMR and MS.
MW	258.1 . 78.2
Purity	≥97% (^{31}P -NMR)
Solubility	Soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)