

# D-myo-Inositol 1,4,5- trisphosphate . hexapotassium salt

Product of phosphatidylinositol phospholipid hydrolysis. Causes release of calcium from intracellular stores.

Citations: 10

[View Online »](#)

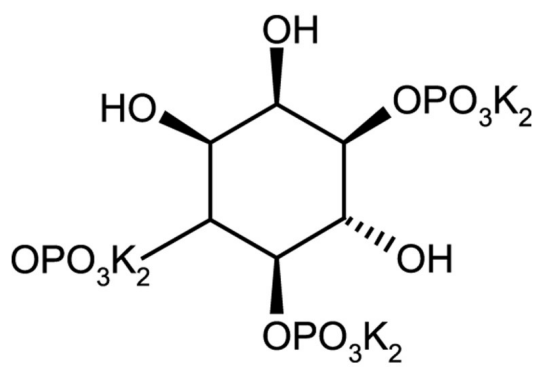
Ordering Information

[Order Online »](#)

ALX-307-007-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Ins(1,4,5)P3 . 6K, IP3
<b>Appearance</b>	White to off-white powder.
<b>CAS</b>	103476-24-0
<b>Formula</b>	$C_6H_9O_{15}P_3 \cdot 6K$
<b>MW</b>	414.1 . 234.6
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	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-  
be@enzolifesciences.com](mailto:info-be@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)