

PRODUCT INFORMATION



D-myo-Inositol-2,3,5,6-tetraphosphate (sodium salt)

Item No. 10008455

Formal Name: D-myo-inositol-2,3,5,6-tetrakis (dihydrogen phosphate), tetrasodium salt

Synonyms: 2,3,5,6-IP₄ (sodium salt), Ins(2,3,5,6)-P₄ (sodium salt)

MF: C₆H₁₂O₁₈P₄ • 4Na

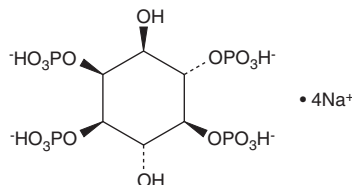
FW: 588.0

Purity: ≥98%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

D-myo-Inositol-2,3,5,6-tetraphosphate (Ins(2,3,5,6)-P₄) (sodium salt) is supplied as a lyophilized powder. Ins(2,3,5,6)-P₄ (sodium salt) is sparingly soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. For biological experiments, we suggest that organic solvent-free aqueous solutions of Ins(2,3,5,6)-P₄ (sodium salt) be prepared by directly dissolving the lyophilized powder in water. The solubility of Ins(2,3,5,6)-P₄ (sodium salt) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ins(2,3,5,6)-P₄ (sodium salt) is produced from Ins(1,2,3,5,6)P₅ via a minor enzymatic phytate degradation pathway by *K. terrigena*.¹ Ins(2,3,5,6)-P₄ (sodium salt) is moderately effective in opening calcium channels relative to Ins(1,4,5)P₃.²

References

1. Greiner, R. and Carlsson, N.-G. *myo*-Inositol phosphate isomers generated by the action of a phytate-degrading enzyme from *Klebsiella terrigena* on phytate. *Can. J. Microbiol.* **52**, 759-768 (2006).
2. DeLisle, S., Radenberg, T., Wintermantel, M.R., *et al.* Second messenger specificity of the inositol trisphosphate receptor: Reappraisal based on novel inositol phosphates. *Am. J. Physiol. Cell Physiol.* **35**, C429-C436 (1994).

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, 07/13/2023

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