

PRODUCT INFORMATION



D-*myo*-Inositol-1,2,3,5-tetraphosphate (sodium salt)

Item No. 10008449

Formal Name: D-*myo*-inositol-1,2,3,5-tetrakis (dihydrogen phosphate), tetrasodium salt

Synonyms: Ins(1,2,3,4)-P₄ (sodium salt), 1,2,3,5-IP₄ (sodium salt)

MF: C₆H₁₂O₁₈P₄ • 4Na

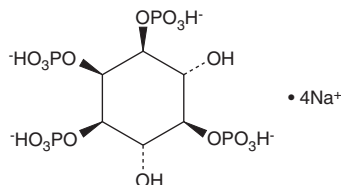
FW: 588.0

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-1,2,3,5-tetraphosphate (Ins(1,2,3,5)-P₄) (sodium salt) is supplied as a lyophilized powder. Ins(1,2,3,5)P₄ (sodium salt) is sparingly soluble in organic solvents. For biological experiments, we suggest that aqueous solutions of Ins(1,2,3,5)-P₄ (sodium salt) be prepared by directly dissolving the lyophilized powder in water. The solubility of Ins(1,2,3,5)-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(1,2,3,5)-P₄ is one of the many inositol phosphate isomers that could act as small, soluble second messengers in the transmission of cellular signals.¹⁻³ Ins(1,2,3,5)-P₄ exhibits relatively weak agonist activity at the Ins(1,4,5)-P₃ receptor expressed in CHO cells, where it induces Ca²⁺ mobilization with an ED₅₀ value of 93.3 μM, making it nearly 1,000-fold less effective than Ins(1,4,5)-P₃ in the same assay.⁴

References

1. Berridge, M.J. Inositol trisphosphate and calcium signalling. *Nature* **361**, 315-325 (1993).
2. Majerus, P.W. Inositol phosphate biochemistry. *Annu. Rev. Biochem.* **61**, 225-250 (1992).
3. Shears, S.B. The versatility of inositol phosphates as cellular signals. *Biochim. Biophys. Acta* **1436**, 49-67 (1998).
4. Burford, N.T., Nahorski, S.R., Chung, S.-K., *et al.* Binding and activity of the nine possible regioisomers of *myo*-inositol tetrakisphosphate at the inositol 1,4,5-trisphosphate receptor. *Cell Calcium* **21**, 301-310 (1997).

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