

PRODUCT INFORMATION

D-*myo*-Inositol-1,2,4,5,6-pentaphosphate (sodium salt)

Item No. 10008452

Formal Name: D-*myo*-inositol-1,2,4,5,6-penta (hydrogen phosphate), pentasodium salt

Synonyms: Ins(1,2,4,5,6)P₅ (sodium salt), 1,2,4,5,6-IP₅ (sodium salt)

MF: C₆H₁₂O₂₁P₅ • 5Na

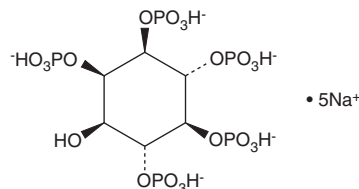
FW: 690.0

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-1,2,4,5,6-pentaphosphate (Ins(1,2,4,5,6)P₅) (sodium salt) is supplied as a lyophilized powder. Ins(1,2,4,5,6)P₅ (sodium salt) is sparingly soluble in organic solvents. For biological experiments, we suggest that aqueous solutions of Ins(1,2,4,5,6)P₅ (sodium salt) be prepared by directly dissolving the lyophilized powder in water. The solubility of Ins(1,2,4,5,6)P₅ (sodium salt) in water is approximately 50 mg/ml.

Description

Ins(1,2,4,5,6)P₅ is one of several different inositol oligophosphate isomers implicated in signal transduction. In human Jurkat T-lymphocytes, as well as in cytosolic extracts from fetal calf thymus, Ins(1,2,4,5,6)P₅ is phosphorylated by a 1/3 kinase to InsP₆.^{1,2}

References

1. Rudolf, M.T., Kaiser, T., Guse, A.H., *et al.* Synthesis and metabolism of the *myo*-inositol pentakisphosphates. *Liebigs Ann.* 1861-1869 (1997).
2. Guse, A.H. and Emmrich, F. T-cell receptor-mediated metabolism of inositol polyphosphates in Jurkat T-lymphocytes. *J. Biol. Chem.* **266**(36), 24498-24502 (1991).

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