

PRODUCT INFORMATION



L-*myo*-Inositol-1,4,5-triphosphate (sodium salt)

Item No. 10008426

Formal Name: 3,5,6-*tris*(dihydrogen phosphate) D-*myo*-inositol, trisodium salt

Synonyms: D-*myo*-Inositol-3,5,6-triphosphate, Ins(1,4,5)P₃, 1,4,5-IP₃

MF: C₆H₁₂O₁₅P₃ • 3Na

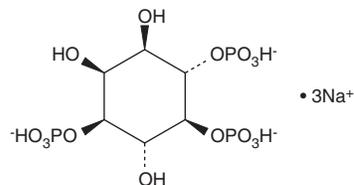
FW: 486.0

Purity: ≥98%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

L-*myo*-Inositol-1,4,5-triphosphate (Ins(1,4,5)P₃) (sodium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ins(1,4,5)P₃ (sodium salt) in water. The solubility of Ins(1,4,5)P₃ (sodium salt) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ins(1,4,5)P₃ is an isomer of the biologically important D-*myo*-inositol-1,4,5-triphosphate. Unlike its isomer, Ins(1,4,5)P₃ does not evoke a rise in intracellular calcium when added to cells.¹ It is not known if Ins(1,4,5)P₃ can act as a competitive inhibitor of biologically-active inositol phosphates.

Reference

1. Denis, G.V. and Ballou, C.E. The Ca²⁺ release activities of D-*myo*-inositol 1,4,5-triphosphate analogs are quantized. *Cell Calcium* **12**(6), 395-401 (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, 12/09/2021

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