

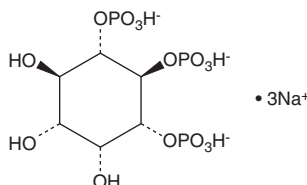
PRODUCT INFORMATION



D-*myo*-Inositol-1,5,6-triphosphate (sodium salt)

Item No. 10008422

CAS Registry No.: 120965-76-6
Formal Name: D-*myo*-inositol-1,5,6-*tris*
(hydrogen phosphate), trisodium salt
Synonyms: Ins(1,5,6)-P₃ (sodium salt), 1,5,6-IP₃
(sodium salt)
MF: C₆H₁₂O₁₅P₃ • 3Na
FW: 486.0
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-1,5,6-triphosphate (Ins(1,5,6)-P₃) (sodium salt) is supplied as a lyophilized powder. For biological experiments, we suggest that organic solvent-free aqueous solutions of Ins(1,5,6)-P₃ (sodium salt) be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of Ins(1,5,6)-P₃ (sodium salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

The inositol phosphates are a family of mono- to poly-phosphorylated compounds that act as messengers, regulating cellular functions including cell cycling, apoptosis, differentiation, and motility.^{1,2} Ins(1,5,6)-P₃ is an intermediate compound, produced by the dephosphorylation of various inositol-tetrakisphosphate forms. The triphosphate can be further metabolized to produce inositol-biphosphate mediators. The biological roles of Ins(1,5,6)-P₃ remain to be determined.

References

1. Majerus, P.W. Inositol phosphate biochemistry. *Annu. Rev. Biochem.* **61**, 225-250 (1992).
2. Berridge, M.J. Inositol trisphosphate and calcium signalling. *Nature* **361**, 315-325 (1993).

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, 07/12/2023

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