

# PRODUCT INFORMATION



## D-chiro-Inositol 1,2,3,4,5,6-hexakisphosphate (sodium salt)

Item No. 9002341

**Formal Name:** (1R,2R,3S,4S,5S,6S)-cyclohexane-(1,2,3,4,5,6)-hexayl hexakis(hydrogen phosphate), hexasodium salt

**MF:**  $C_6H_{12}O_{24}P_6 \cdot 6Na$

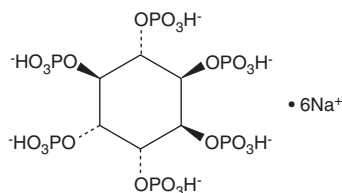
**FW:** 791.9

**Purity:**  $\geq 98\%$

**Supplied as:** A crystalline solid

**Storage:**  $-20^{\circ}C$

**Stability:**  $\geq 5$  years



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

### Laboratory Procedures

D-chiro-Inositol 1,2,3,4,5,6-hexakisphosphate (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the D-chiro-inositol 1,2,3,4,5,6-hexakisphosphate (sodium salt) in water. The solubility of D-chiro-inositol 1,2,3,4,5,6-hexakisphosphate (sodium salt) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

D-chiro-Inositol 1,2,3,4,5,6-hexakisphosphate is a stereoisomer of inositol hexakisphosphate ( $IP_6$ ), an organic phosphorous compound found in the environment.<sup>1</sup> D-chiro-Inositol 1,2,3,4,5,6-hexakisphosphate has been identified in soil samples from Madagascar, New Zealand, and the Falkland Islands and in freshwater sediment from Danish lakes.<sup>2,3</sup>

### References

1. Turner, B.L., Cheesman, A.W., Godage, H.Y., *et al.* Determination of *neo*- and D-chiro-inositol hexakisphosphate in soils by solution  $^{31}P$  NMR spectroscopy. *Environ. Sci. Technol.* **46**(9), 4994-5002 (2012).
2. Jørgensen, C., Jensen, H.S., Andersen, F.Ø., *et al.* Occurrence of orthophosphate monoesters in lake sediments: Significance of *myo*- and *scyllo*-inositol hexakisphosphate. *J. Environ. Monit.* **13**(8), 2328-2334 (2011).
3. Jørgensen, C., Turner, B.L., and Reitzel, K. Identification of inositol hexakisphosphate binding sites in soils by selective extraction and solution  $^{31}P$  NMR spectroscopy. *Geoderma* **257-258**, 22-28 (2015).

#### 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, 07/25/23

#### 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](http://WWW.CAYMANCHEM.COM)