

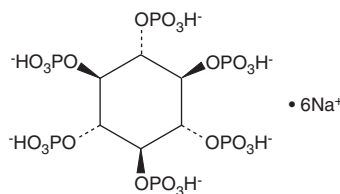
# PRODUCT INFORMATION



## scyllo-Inositol hexakisphosphate (sodium salt)

Item No. 9002339

**Formal Name:** scyllo-inositol, 1,2,3,4,5,6-hexakis(dihydrogen phosphate), hexasodium salt  
**Synonym:** scyllo-IP<sub>6</sub>  
**MF:** C<sub>6</sub>H<sub>12</sub>O<sub>24</sub>P<sub>6</sub> • 6Na  
**FW:** 791.9  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Laboratory Procedures

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

### Description

scyllo-Inositol hexakisphosphate is a stereoisomer of inositol hexakisphosphate (IP<sub>6</sub>), an organic phosphorous compound found in the environment.<sup>1</sup> scyllo-Inositol 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> scyllo-Inositol hexakisphosphate has been used as a negative control for the stereospecific effects of myo-inositol hexakisphosphate on I<sub>K,in</sub> in intact guard cells.<sup>4</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 <sup>31</sup>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 <sup>31</sup>P NMR spectroscopy. *Geoderma* **257-258**, 22-28 (2015).
4. Lemtiri-Chlieh, F., MacRobbie, E.A., and Brearley, C.A. Inositol hexakisphosphate is a physiological signal regulating the K<sup>+</sup>-inward rectifying conductance in guard cells. *Proc. Natl. Acad. Sci. U.S.A.* **97(15)**, 8687-8692 (2000).

#### 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, 07/24/2023

#### 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