

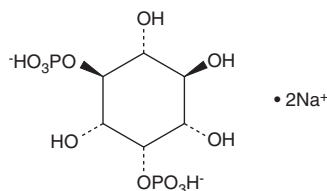
# PRODUCT INFORMATION



## D-*myo*-Inositol-2,4-diphosphate (sodium salt)

Item No. 10008419

**Formal Name:** D-*myo*-inositol-2,4-*bis*(hydrogen phosphate), disodium salt  
**Synonyms:** Ins(2,4)-P<sub>2</sub> (sodium salt), 2,4-IP<sub>2</sub> (sodium salt)  
**MF:** C<sub>6</sub>H<sub>12</sub>O<sub>12</sub>P<sub>2</sub> • 2Na  
**FW:** 384.1  
**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-2,4-diphosphate (Ins(2,4)-P<sub>2</sub>) (sodium salt) is supplied as a lyophilized powder. For biological experiments, we suggest that organic solvent-free aqueous solutions of Ins(2,4)-P<sub>2</sub> (sodium salt) be prepared by directly dissolving the lyophilized powder in aqueous buffers. The solubility of Ins(2,4)-P<sub>2</sub> (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.<sup>1,2</sup> Ins(2,4)-P<sub>2</sub> is an intermediate compound, produced by the dephosphorylation of either inositol-2,3,4-triphosphate or inositol-2,4,5-triphosphate. The diphosphate can be further metabolized to produce inositol-2-monophosphate. The biological roles of Ins(2,4)-P<sub>2</sub> 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