

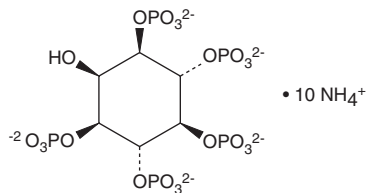
PRODUCT INFORMATION



D-*myo*-Inositol-1,3,4,5,6-pentaphosphate (ammonium salt)

Item No. 10009851

Formal Name:	D- <i>myo</i> -inositol-1,3,4,5,6-penta (hydrogen phosphate), decaammonium salt
Synonyms:	Ins(1,3,4,5,6)P ₅ , 1,3,4,5,6-IP ₅ (sodium salt)
MF:	C ₆ H ₇ O ₂₁ P ₅ • 10NH ₄
FW:	750.4
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-1,3,4,5,6-pentaphosphate (Ins(1,3,4,5,6)P₅) (ammonium salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the Ins(1,3,4,5,6)P₅ (ammonium salt) in water. The solubility of Ins(1,3,4,5,6)P₅ (ammonium salt) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Ins(1,3,4,5,6)P₅ is one of the many inositol phosphate isomers that act as small, soluble second messengers in the transmission of cellular signals.¹⁻³ It can be interconverted with Ins(3,4,5,6)P₄ by a 1-kinase/1-phosphatase cycle, as well as with Ins(1,4,5,6)P₄ in a 3-kinase/3-phosphatase cycle.³ Ins(1,3,4,5,6)P₅ inhibits the phosphorylation and kinase activity of Akt/PKB, inducing apoptosis in ovarian, lung, and breast cancer cells.⁴ It exhibits antiangiogenic activity *in vitro*, blocking capillary tube formation of HUVEC, as well as antitumor effects against cancer xenografts in nude mice.⁵ Ins(1,3,4,5,6)P₅ binds to the PH domain of Grp1 with a K_d of 590 nM.⁶

References

1. Berridge, M.J. Inositol trisphosphate and calcium signalling. *Nature* **361(6410)**, 315-325 (1993).
2. Majerus, P.W. Inositol phosphate biochemistry. *Annu. Rev. Biochem.* **61**, 225-250 (1992).
3. Shears, S.B. The versatility of inositol phosphates as cellular signals. *Biochim. Biophys. Acta.* **1436(1-2)**, 49-67 (1998).
4. Piccolo, E., Vignati, S., Maffucci, T., *et al.* Inositol pentakisphosphate promotes apoptosis through the PI 3-K/Akt pathway. *Oncogene* **23(9)**, 1754-1765 (2004).
5. Maffucci, T., Piccolo, E., Cumashi, A., *et al.* Inhibition of the phosphatidylinositol 3-kinase/Akt pathway by inositol pentakisphosphate results in antiangiogenic and antitumor effects. *Cancer Res.* **65(18)**, 8339-8349 (2005).
6. Kavran, J.M., Klein, D.E., Lee, A., *et al.* Specificity and promiscuity in phosphoinositide binding by pleckstrin homology domains. *J. Biol. Chem.* **273(46)**, 30497-30508 (1998).

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, 07/13/2023

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