

PRODUCT INFORMATION



PTP1B-IN-2

Item No. 36824

CAS Registry No.: 1919853-46-5

Formal Name: 5-[[[4-[(2-methoxy-2-oxoethyl)(methylsulfonyl)amino]phenyl]methyl][(phenylmethyl)sulfonyl]amino]-2-[(4-methylphenyl)methoxy]-benzoic acid, methyl ester

Synonyms: Protein Tyrosine Phosphatase 1B Inhibitor 2, PTPN1 Inhibitor 2, Tyrosine-protein Phosphatase Non-Receptor Type 1 Inhibitor 2

MF: $C_{34}H_{36}N_2O_9S_2$

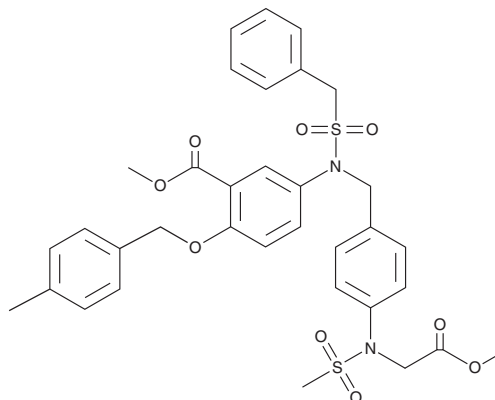
FW: 680.8

Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years



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

Laboratory Procedures

PTP1B-IN-2 is supplied as a solid. A stock solution may be made by dissolving the PTP1B-IN-2 in the solvent of choice, which should be purged with an inert gas. PTP1B-IN-2 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PTP1B-IN-2 in these solvents is approximately 10 mg/ml.

Description

PTP1B-IN-2 is an inhibitor of protein tyrosine phosphatase 1B (PTP1B; $IC_{50} = 50$ nM).¹ It is selective for PTP1B over T cell protein tyrosine phosphatase (TCPTP), Src homology region 2 domain-containing phosphatase 1 (SHP-1), SHP-2, and leukocyte common antigen-related (LAR) phosphatase (IC_{50} s = 760, 2,450, 2,160, and 12,610 nM, respectively). PTP1B-IN-2 (5-20 μM) enhances insulin-induced glucose uptake in L6 myotubes.

Reference

1. Liu, P., Du, Y., Song, L., *et al.* Discovery of novel, high potent, ABC type PTP1B inhibitors with TCPTP selectivity and cellular activity. *Eur. J. Med. Chem.* **118**, 27-33 (2016).

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, 11/01/2023

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