

PRODUCT INFORMATION



PTP Inhibitor IV

Item No. 18458

CAS Registry No.: 329317-98-8

Formal Name: N,N'-[1,4-phenylenebis[(1-methylethylidene)-4,1-phenylene]]bis[1,1,1-trifluoromethanesulfonamide

Synonym: Protein Tyrosine Phosphatase Inhibitor IV

MF: C₂₆H₂₆F₆N₂O₄S₂

FW: 608.6

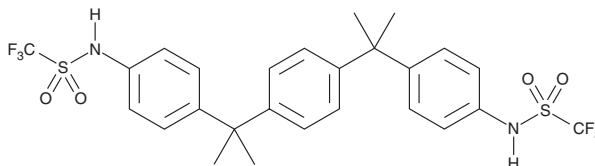
Purity: ≥98%

UV/Vis.: λ_{max}: λ_{max}: 229 nm

Supplied as: A crystalline 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

PTP inhibitor IV is supplied as a crystalline solid. A stock solution may be made by dissolving the PTP inhibitor IV in the solvent of choice, which should be purged with an inert gas. PTP inhibitor IV is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PTP inhibitor IV in ethanol and DMSO is approximately 25 mg/ml and approximately 30 mg/ml in DMF.

PTP inhibitor IV is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PTP inhibitor IV should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PTP inhibitor IV has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Protein tyrosine phosphatases (PTPs) remove phosphate from tyrosine residues of cellular proteins. Reversible phosphorylation catalyzed by the coordinated actions of protein tyrosine kinases and phosphatases is key to the regulation of the signaling events that control cell growth and proliferation, differentiation, and survival or apoptosis, as well as adhesion and motility. PTP inhibitor IV is an uncharged, 1,4-di-substituted, phenyl-linked bis-trifluoromethylsulfonamido phosphate mimetic that acts as a reversible, competitive, and active-site directed inhibitor of SHP-2, PTP1B, PTP-ε, PTP-Meg-2, PTP-σ, PTP-β, and PTP-μ (IC₅₀s = 1.8, 2.5, 8.4, 13, 20, 6.4, and 6.7 μM, respectively).¹

Reference

- Huang, P., Ramphal, J., Wei, J., *et al.* Structure-based design and discovery of novel inhibitors of protein tyrosine phosphatases. *Bioor. Med. Chem.* **11**, 1835-1849 (2003).

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, 10/13/2022

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