

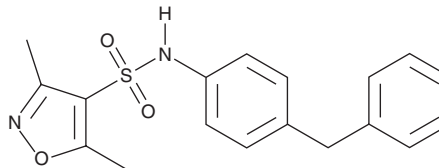
PRODUCT INFORMATION



PP5-IN-1

Item No. 43071

CAS Registry No.: 1022417-69-1
Formal Name: 3,5-dimethyl-N-[4-(phenylmethyl)phenyl]-4-isoxazolesulfonamide
Synonyms: P053, PP5-Inhibitor 1, Protein Phosphatase 5 Inhibitor 1
MF: C₁₈H₁₈N₂O₃S
FW: 342.4
Purity: ≥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

PP5-IN-1 is supplied as a solid. A stock solution may be made by dissolving the PP5-IN-1 in the solvent of choice, which should be purged with an inert gas. PP5-IN-1 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

PP5-IN-1 is an inhibitor of protein phosphatase 5 (PP5; IC₅₀ = 0.706 μM).¹ It induces apoptosis in 786-O renal cell carcinoma (RCC) cells when used at a concentration of 10 μM.

Reference

1. Ahanin, E.F., Sager, R.A., Backe, S.J., et al. Catalytic inhibitor of Protein Phosphatase 5 activates the extrinsic apoptotic pathway by disrupting complex II in kidney cancer. *Cell Chem. Biol.* **30(10)**, 1223-1234 (2023).

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, 04/09/2025

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