

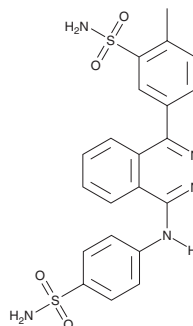
PRODUCT INFORMATION



ENPP3 Inhibitor

Item No. 37760

CAS Registry No.: 333413-02-8
Formal Name: 5-[4-[[4-(aminosulfonyl)phenyl]amino]-1-phthalazinyl]-2-methylbenzenesulfonamide
MF: C₂₁H₁₉N₅O₄S₂
FW: 469.5
Purity: ≥95%
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

ENPP3 inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the ENPP3 inhibitor in the solvent of choice, which should be purged with an inert gas. ENPP3 inhibitor is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

Description

ENPP3 inhibitor is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase family member 3 (ENPP3; K_i = 53.7 nM).¹ It is selective for ENPP3 over ENPP1, ENPP2, tissue-nonspecific alkaline phosphatase (TNAP), ecto-5'-nucleotidase (CD73), nucleoside triphosphate diphosphohydrolase 1 (NTPDase 1), NTPDase 2, and NTPDase 3. ENPP3 inhibitor also inhibits carbonic anhydrase II (CAII) and CAIX (K_is = 74.7 and 20.3 nM, respectively).

Reference

1. Lee, S.-Y., Namasivayam, V., Boshta, N.M., *et al.* Discovery of potent nucleotide pyrophosphatase/phosphodiesterase3 (NPP3) inhibitors with ancillary carbonic anhydrase inhibition for cancer (immuno) therapy. *RSC Med. Chem.* **12**(7), 1187-1206 (2021).

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, 10/09/2025

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