

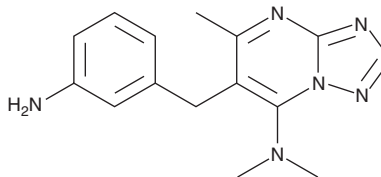
PRODUCT INFORMATION



ENPP1 Inhibitor C

Item No. 29809

CAS Registry No.: 2378640-92-5
Formal Name: 6-[(3-aminophenyl)methyl]-N,N,5-trimethyl-[1,2,4]triazolo[1,5-a]pyrimidin-7-amine
MF: C₁₅H₁₈N₆
FW: 282.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

ENPP1 inhibitor C is supplied as a crystalline solid. A stock solution may be made by dissolving the ENPP1 inhibitor C in the solvent of choice, which should be purged with an inert gas. ENPP1 inhibitor C is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of ENPP1 inhibitor C in these solvents is approximately 30 mg/ml.

Description

ENPP1 inhibitor C is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase 1 (ENPP1; IC₅₀ = 0.26 μM in a cell-free assay).¹ It is selective for ENPP1 over ENPP2-7 at 10 μM. ENPP1 inhibitor C decreases ENPP1 activity in MDA-MB-231 human breast and C6 rat glioma cancer cells when used at a concentration of 10 μM.

Reference

1. Kawaguchi, M., Han, X., Hisada, T., *et al.* Development of an ENPP1 fluorescence probe for inhibitor screening, cellular imaging, and prognostic assessment of malignant breast cancer. *J. Med. Chem.* **62(20)**, 9254-9269 (2019).

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, 02/13/2020

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