

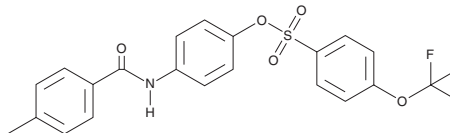
PRODUCT INFORMATION



ENPP3 Inhibitor 4g

Item No. 37647

CAS Registry No.: 2803505-78-2
Formal Name: 4-(trifluoromethoxy)-benzenesulfonic acid, 4-[(4-methylbenzoyl)amino]phenyl ester
MF: C₂₁H₁₆F₃NO₅S
FW: 451.4
Purity: ≥95%
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

ENPP3 inhibitor 4g is supplied as a solid. A stock solution may be made by dissolving the ENPP3 inhibitor 4g in the solvent of choice, which should be purged with an inert gas. ENPP3 inhibitor 4g is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ENPP3 inhibitor 4g in these solvents is approximately 12 and 16 mg/ml, respectively. ENPP3 inhibitor 4g is slightly soluble in ethanol.

Description

ENPP3 inhibitor 4g is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase family member 1 (ENPP1) and ENPP3 (IC₅₀s = 0.45 and 0.19 μM, respectively).¹

Reference

1. Ullah, S., Pelletier, J., Sévigny, J., *et al.* Synthesis and biological evaluation of arylamide sulphonate derivatives as ectonucleotide pyrophosphatase/phosphodiesterase-1 and -3 inhibitors. *ACS Omega* **7**(30), 26905-26918 (2022).

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, 05/08/2024

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