

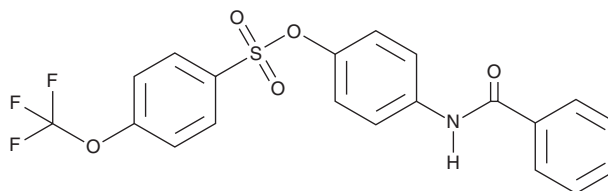
PRODUCT INFORMATION



ENPP3 Inhibitor 4t

Item No. 37646

CAS Registry No.: 2803505-90-8
Formal Name: 4-benzamidophenyl 4-(trifluoromethoxy) benzenesulfonate
MF: C₂₀H₁₄F₃NO₅S
FW: 437.4
Purity: ≥95%
UV/Vis.: λ_{max}: 270 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

ENPP3 inhibitor 4t is supplied as a crystalline solid. A stock solution may be made by dissolving the ENPP3 inhibitor 4t in the solvent of choice, which should be purged with an inert gas. ENPP3 inhibitor 4t is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ENPP3 inhibitor 4t in these solvents is approximately 3 mg/ml.

Description

ENPP3 inhibitor 4t is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase family member 3 (ENPP3; IC₅₀ = 0.15 μM).¹ It is selective for ENPP3 over ENPP1 (IC₅₀ = 41.4 μM).

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, 09/04/2024

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