

PRODUCT INFORMATION



ENPP3 Inhibitor 5J

Item No. 40265

Formal Name: 2-(2-(1H-indol-3-yl)acetamido)phenyl trifluoromethanesulfonate

MF: C₁₇H₁₃F₃N₂O₄S

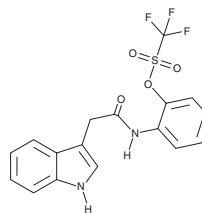
FW: 398.4

Purity: ≥95%

Supplied as: A 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

ENPP3 inhibitor 5J is supplied as a solid. A stock solution may be made by dissolving the ENPP3 inhibitor 5J in the solvent of choice, which should be purged with an inert gas. ENPP3 inhibitor 5J is soluble in acetonitrile and chloroform.

Description

ENPP3 inhibitor 5J is an inhibitor of ectonucleotide pyrophosphatase/phosphodiesterase family member 1 (ENPP1) and ENPP3 (IC₅₀s = 0.59 and 0.62 μM, respectively, for the human enzymes).¹

Reference

1. Khan Jadoon, M.S., Pelletier, J., Sévigny, J., *et al.* Synthesis of new class of indole acetic acid sulfonate derivatives as ectonucleotidases inhibitors. *RSC Adv.* **13(42)**, 29496-29511 (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, 02/08/2024

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