

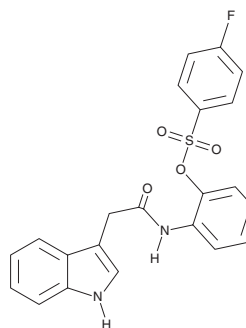
PRODUCT INFORMATION



h-e5'NT Inhibitor 5C

Item No. 40266

Formal Name: 2-(2-(1H-indol-3-yl)acetamido)phenyl
4-fluorobenzenesulfonate
Synonym: Human Ecto-5'-nucleotidase Inhibitor 5C
MF: C₂₂H₁₇FN₂O₄S
FW: 424.4
Purity: ≥95%
Supplied as: A neat 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

h-e5'NT Inhibitor 5C is supplied as a neat solid. A stock solution may be made by dissolving the *h*-e5'NT inhibitor 5C in the solvent of choice, which should be purged with an inert gas. *h*-e5'NT Inhibitor 5C is soluble in acetonitrile, methanol, and DMSO.

Description

h-e5'NT Inhibitor 5C is an inhibitor of ecto-5'-nucleotidase, also known as CD73 (IC₅₀s = 0.37 and 1.66 μM for the human and rat enzymes, respectively).¹ It also inhibits ectonucleotide pyrophosphate/phosphodiesterase family member 1 (ENPP1), ENPP3, and tissue-non-specific alkaline phosphatase (TNAP; IC₅₀s = 1.12, 1.37, and 1.48 μM, respectively, for the human enzymes).

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, 02/06/2025

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