

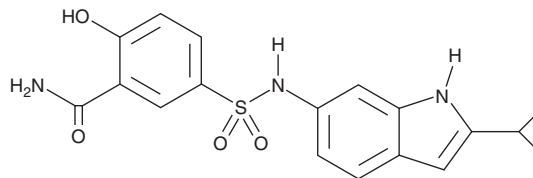
PRODUCT INFORMATION



CD73-IN-1

Item No. 30959

CAS Registry No.: 2132396-40-6
Formal Name: 5-[[[(2-cyclopropyl-1H-indol-6-yl)amino]sulfonyl]-2-hydroxy-benzamide
MF: C₁₈H₁₇N₃O₄S
FW: 371.4
Purity: ≥98%
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

CD73-IN-1 is supplied as a solid. A stock solution may be made by dissolving the CD73-IN-1 in the solvent of choice, which should be purged with an inert gas. CD73-IN-1 is soluble in DMSO.

Description

CD73-IN-1 is an inhibitor of ecto-5'-nucleotidase, also known as CD73, with an IC₅₀ value of ≤316.23 nM for the human enzyme in a cell-free assay.¹

Reference

1. Adams, J.L., Duffy, K.J. 5-sulfamoyl-2-hydroxybenzamide derivatives. *GlaxoSmithKline*. **WO2017/153952A1** (2017).

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, 11/04/2022

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