

PRODUCT INFORMATION

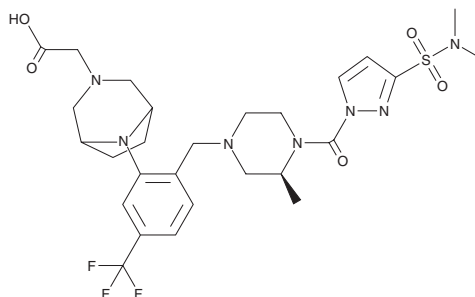


ABD957

Item No. 44290

CAS Registry No.: 3007772-60-0
Formal Name: 8-[2-[[[(3S)-4-[[3-[(dimethylamino)sulfonyl]-1H-pyrazol-1-yl]carbonyl]-3-methyl-1-piperazinyl]methyl]-5-(trifluoromethyl)phenyl]-3,8-diazabicyclo[3.2.1]octane-3-acetic acid

MF: C₂₇H₃₆F₃N₇O₅S
FW: 627.7
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

ABD957 is supplied as a solid. A stock solution may be made by dissolving the ABD957 in the solvent of choice, which should be purged with an inert gas. ABD957 is soluble in DMSO.

Description

ABD957 is an inhibitor of α/β -hydrolase domain-containing protein 17 (ABHD17; IC₅₀ = 0.21 μ M for ABHD17B).¹ It is selective for ABHD17A-C over a panel of 52 serine hydrolases, including the depalmitoylases ABHD10, palmitoyl-protein thioesterase 1 (PPT1), and lysophospholipase 1 (LYPLA1) and LYPLA2, at 10 μ M but does inhibit ABHD6, ABHD11, ABHD13, and carboxylesterase 1 (CES1) and CES2 at the same concentration. ABD957 increases the palmitoylation of N-Ras in OCI-AML-3 acute myeloid leukemia (AML) cells expressing GFP-N-Ras^{G12D} (EC₅₀ = 29 nM). It also increases the plasma membrane localization of N-Ras in the same cells when used at a concentration of 500 nM. ABD957 inhibits the growth of OCI-AML-3 cells in a concentration-dependent manner.

Reference

1. Remsberg, J.R., Suciu, R.M., Zambetti, N.A., *et al.* ABHD17 regulation of plasma membrane palmitoylation and N-Ras-dependent cancer growth. *Nat. Chem. Biol.* **17**(8), 856-864 (2021).

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, 09/23/2025

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