

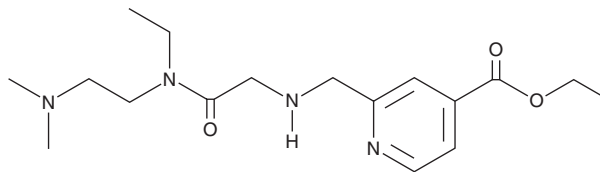
PRODUCT INFORMATION



KDM5-C70

Item No. 36817

CAS Registry No.: 1596348-32-1
Formal Name: 2-[[[2-[[2-(dimethylamino)ethyl]ethylamino]-2-oxoethyl]amino]methyl]-4-pyridinecarboxylic acid, ethyl ester
Synonym: KDOAM-21
MF: C₁₇H₂₈N₄O₃
FW: 336.4
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

KDM5-C70 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used.

Description

KDM5-C70 is an inhibitor of Jumonji AT rich interactive domain 1B (JARID1B), also known as lysine-specific demethylase 5B (KDM5B; IC₅₀ = 49 nM).¹ It is selective for JARID1B over KDM4C (IC₅₀ = 2,400 nM).

Reference

1. Tumber, A., Nuzzi, A., Hookway, E.S., *et al.* Potent and selective KDM5 inhibitor stops cellular demethylation of H3K4me3 at transcription start sites and proliferation of MM1S myeloma cells. *Cell Chem. Biol.* **24**(3), 371-380 (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, 12/21/2022

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