

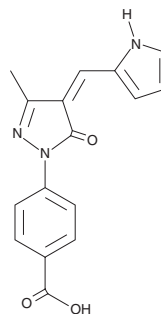
PRODUCT INFORMATION



KAT8 Inhibitor 19

Item No. 38977

CAS Registry No.: 28532-21-0
Formal Name: 4-[4,5-dihydro-3-methyl-5-oxo-4-(1H-pyrrol-2-ylmethylene)-1H-pyrazol-1-yl]-benzoic acid
Synonym: Lysine Acetyltransferase 8 Inhibitor 19
MF: C₁₆H₁₃N₃O₃
FW: 295.3
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

KAT8 inhibitor 19 is supplied as a solid. A stock solution may be made by dissolving the KAT8 inhibitor 19 in the solvent of choice, which should be purged with an inert gas. KAT8 inhibitor 19 is soluble in acetonitrile and DMSO.

Description

KAT8 inhibitor 19 is an inhibitor of lysine acetyltransferase 8 (KAT8; IC₅₀ = 12.1 μM).¹ It is selective for KAT8 over KAT2A, KAT2B, KAT3B, KAT6A, and KAT7 at 200 μM. KAT8 (50 μM) reduces histone 4 lysine 16 acetylation (H4K16Ac) in HT-29 colorectal carcinoma cells. It reduces the viability of HCT116, H1299, A549, and U937 cancer cells (IC₅₀s = 39.4, 52.1, 41, and 30.1 μM, respectively).

Reference

1. Fiorentino, F., Sementilli, S., Menna, M., *et al.* First-in-class selective inhibitors of the lysine acetyltransferase KAT8. *J. Med. Chem.* **66**(10), 6591-6616 (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, 03/05/2024

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