

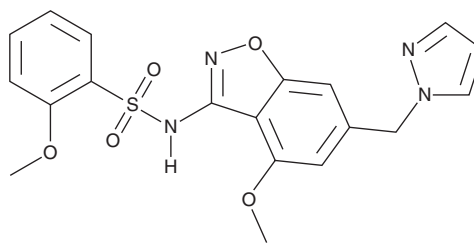
PRODUCT INFORMATION



PF-07248144

Item No. 43955

CAS Registry No.: 2569008-99-5
Formal Name: 2-methoxy-N-[4-methoxy-6-(1H-pyrazol-1-ylmethyl)-1,2-benzisoxazol-3-yl]-benzenesulfonamide
MF: C₁₉H₁₈N₄O₅S
FW: 414.4
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

PF-07248144 is supplied as a solid. A stock solution may be made by dissolving the PF-07248144 in the solvent of choice, which should be purged with an inert gas. PF-07248144 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

PF-07248144 is an inhibitor of lysine acetyltransferase 6A (KAT6A), also known as monocytic leukemia zinc finger protein (MOZ) and KAT6B, also known as monocytic leukemia zinc finger protein-related factor (MORF).¹

Reference

1. Mukohara, T., Park, Y.H., Sommerhalder, D., et al. Inhibition of lysine acetyltransferase KAT6 in ER⁺HER2⁻ metastatic breast cancer: A phase 1 trial. *Nat. Med.* **30(8)**, 2242-2250 (2024).

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, 01/16/2025

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