

PRODUCT INFORMATION



FHT-2344

Item No. 42043

CAS Registry No.: 2468058-90-2

Formal Name: N-[(1S)-1-(methoxymethyl)-2-[[4-[3-(1-methyl-1H-pyrazol-3-yl)phenyl]-2-thiazolyl]amino]-2-oxoethyl]-1-(methylsulfonyl)-1H-pyrrole-3-carboxamide

MF: C₂₃H₂₄N₆O₅S₂

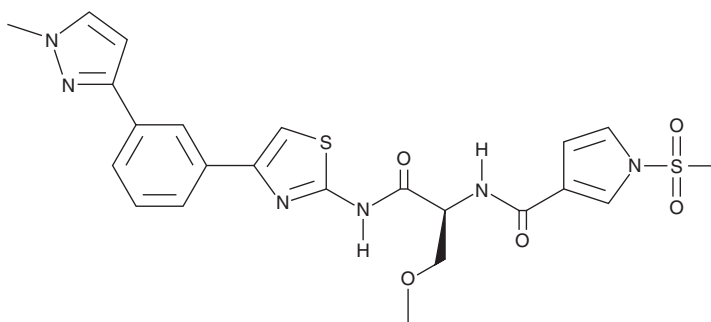
FW: 528.6

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

FHT-2344 is supplied as a solid. A stock solution may be made by dissolving the FHT-2344 in the solvent of choice, which should be purged with an inert gas. FHT-2344 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

FHT-2344 is an inhibitor of Brahma homolog (BRM), also known as SWI/SNF-related, matrix-associated, actin-dependent regulator of chromatin, subfamily A, member 2 (SMARCA2), and Brahma-related gene 1 (BRG1), also known as SMARCA4 (IC₅₀s = ≥0.01 μM for both).¹

Reference

1. Vaswani, R.G., Anthony, N.G., Millan, D. S., *et al.* Compounds and uses thereof. *Foghorn Therapeutics Inc.* WO2020/160180A1 (2020).

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, 10/21/2024

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