

PRODUCT INFORMATION



BI-3231

Item No. 41147

CAS Registry No.: 2894848-07-6

Formal Name: 1-[[5-(2,4-difluoro-3-hydroxyphenyl)-1,3,4-thiadiazol-2-yl]methyl]-3-ethyl-5-methyl-2,4(1H,3H)-pyrimidinedione

MF: C₁₆H₁₄F₂N₄O₃S

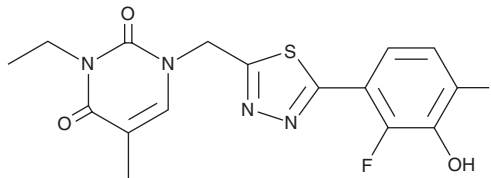
FW: 380.4

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

BI-3231 is supplied as a solid. A stock solution may be made by dissolving the BI-3231 in the solvent of choice, which should be purged with an inert gas. BI-3231 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

BI-3231 is an inhibitor of 17 β -hydroxysteroid dehydrogenase 13 (17 β -HSD13; K_i = 0.7 nM), also known as HSD17B13.¹

Reference

1. Thamm, S., Willwacher, M.K., Aspnes, G.E., *et al.* Discovery of a novel potent and selective HSD17B13 inhibitor, BI-3231, a well-characterized chemical probe available for open science. *J. Med. Chem.* **66**(4), 2832-2850 (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, 01/17/2025

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