

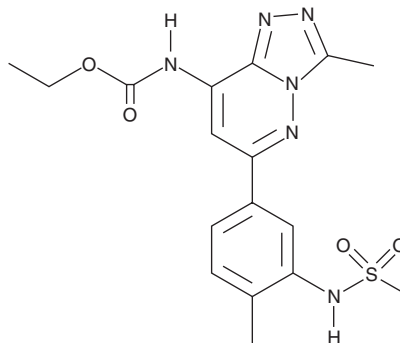
PRODUCT INFORMATION



Bromosporine

Item No. 14119

CAS Registry No.: 1619994-69-2
Formal Name: ethyl (3-methyl-6-(4-methyl-3-(methylsulfonamido)phenyl)-[1,2,4]triazolo[4,3-b]pyridazin-8-yl)carbamate
MF: C₁₇H₂₀N₆O₄S
FW: 404.4
Purity: ≥98%
UV/Vis.: λ_{max}: 255, 293 nm
Supplied as: A crystalline 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

Bromosporine is supplied as a crystalline solid. A stock solution may be made by dissolving the bromosporine in the solvent of choice, which should be purged with an inert gas. Bromosporine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of bromosporine in ethanol is approximately 0.25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Bromosporine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, bromosporine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Bromosporine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Bromosporine is a non-selective bromodomain inhibitor. At 1 μM, it has been shown to accelerate fluorescence recovery after photobleaching (FRAP) recovery of BRD4 and CREBBP bromodomains in HeLa cells.¹ This compound may be used in determining the biological roles of reader domains as well as a tool for the validation of functional assays. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Gallenkamp, D., Gelato, K.A., Haendler, B., *et al.* Bromodomains and their pharmacological inhibitors. *Chem. Med. Chem.* **9**(3), 438-464 (2014).

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, 12/08/2022

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