

PRODUCT INFORMATION



OF-1

Item No. 17124

CAS Registry No.: 919973-83-4
Formal Name: 4-bromo-N-(2,3-dihydro-6-methoxy-1,3-dimethyl-2-oxo-1H-benzimidazol-5-yl)-2-methylbenzenesulfonamide

MF: C₁₇H₁₈BrN₃O₄S

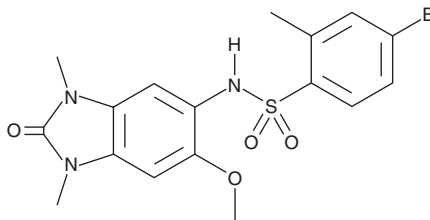
FW: 440.3

Purity: ≥98%

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

OF-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the OF-1 in the solvent of choice, which should be purged with an inert gas. OF-1 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of OF-1 in these solvents is approximately 1 and 5 mg/ml, respectively.

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

Description

OF-1 is a chemical probe for the bromodomains of the Bromodomain and PHD Finger-containing (BRPF) family of proteins. It binds BRPF1B, BRPF2, and BRPF3 with K_d values of 0.1, 0.5, and 2.4 μ M, respectively, as determined by isothermal titration calorimetry. OF-1 increases thermal stability of full length BRPF1B at 1 μ M, in the cellular thermal shift assay. It also demonstrates accelerated fluorescence recovery after photobleaching (FRAP) at 5 μ M in the BRPF2 FRAP assay. OF-1 shows modest general cytotoxicity. See the Structural Genomics Consortium (SGC) website for further information.

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, 11/08/2022

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