

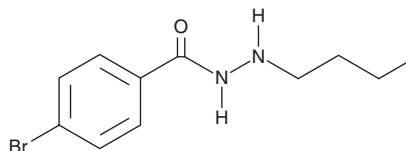
PRODUCT INFORMATION



UF010

Item No. 21273

CAS Registry No.: 537672-41-6
Formal Name: 4-bromo-benzoic acid, 2-butylhydrazide
MF: C₁₁H₁₅BrN₂O
FW: 271.2
Purity: ≥98%
UV/Vis.: λ_{max}: 240 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

UF010 is supplied as a crystalline solid. A stock solution may be made by dissolving the UF010 in the solvent of choice. UF010 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UF010 in these solvents is approximately 30 mg/ml.

UF010 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UF010 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. UF010 has a solubility of approximately 0.12 mg/ml in a 1:7 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UF010 is a class I HDAC inhibitor with IC₅₀ values of 0.06, 0.1, 0.5, and 1.5 μM for HDAC3, 2, 1, and 8, respectively.¹ It is more than 6-fold selective for these class I HDACs over any additional HDACs.¹ UF010 is reported to inhibit cancer cell proliferation with a mean GI₅₀ value of 2.94 μM when screened against the NCI-60 panel of cancer cell lines and functions by blocking the G₁/S phase of the cell cycle.¹

Reference

1. Wang, Y., Stowe, R.L., Pinello, C.E., *et al.* Identification of histone deacetylase inhibitors with benzoylhydrazide scaffold that selectively inhibit class I histone deacetylases. *Chem. Biol.* **22(2)**, 273-284 (2015).

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, 12/22/2022

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