

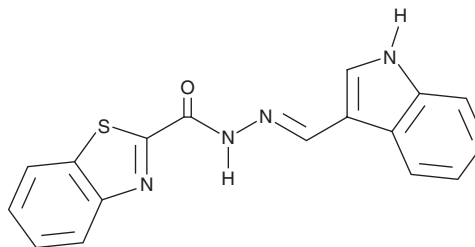
PRODUCT INFORMATION



CAY10777

Item No. 32712

CAS Registry No.: 2460648-40-0
Formal Name: 2-benzothiazolecarboxylic acid, (2E)-2-(1H-indol-3-ylmethylene)hydrazide
MF: C₁₇H₁₂N₄OS
FW: 320.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 260, 358 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10777 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10777 in the solvent of choice, which should be purged with an inert gas. CAY10777 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10777 in these solvents is approximately 10 mg/ml.

CAY10777 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10777 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10777 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10777 is an antiproliferative agent.¹ It inhibits the proliferation of HepG2 cells *in vitro* (IC₅₀ = 0.78 μM).

Reference

1. Liu, K., Ding, Y., and Kang, C. Synthesis and antiproliferative activity of new N-acylhydrazone derivatives containing benzothiazole and indole based moiety. *Pharm. Chem. J.* **54**, 345-352 (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, 11/24/2020

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