

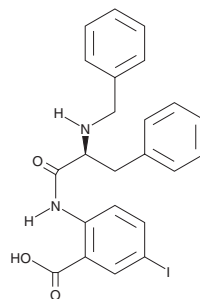
PRODUCT INFORMATION



CW069

Item No. 22936

CAS Registry No.: 1594094-64-0
Formal Name: 5-iodo-2-[[[(2S)-1-oxo-3-phenyl-2-[(phenylmethyl)amino]propyl]amino]-benzoic acid
MF: C₂₃H₂₁IN₂O₃
FW: 500.3
Purity: ≥95%
UV/Vis.: λ_{max}: 265 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

CW069 is supplied as a crystalline solid. A stock solution may be made by dissolving the CW069 in the solvent of choice. CW069 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CW069 in ethanol is approximately 2.5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CW069 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CW069 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CW069 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

CW069 is an allosteric inhibitor of KIFC/HSET (IC₅₀ = 75 μM), a microtubule motor protein involved in bipolar spindle assembly and centrosome clustering.¹ It increases the number of multipolar spindles in N1E-115 mouse neuroblastoma cells with supernumerary centrosomes when used at concentrations of 100 and 200 μM, without affecting the morphology of mitotic spindles in normal human dermal fibroblast (NHDF) cells. CW069 inhibits proliferation of N1E-115, but not NHDF, cells (IC₅₀s = 86 and 181 μM, respectively).

Reference

1. Watts, C.A., Richards, F.M., Bender, A., *et al.* Design, synthesis, and biological evaluation of an allosteric inhibitor of HSET that targets cancer cells with supernumerary centrosomes. *Chem. Biol.* **20(11)**, 1399-1410 (2013).

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, 12/12/2022

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