

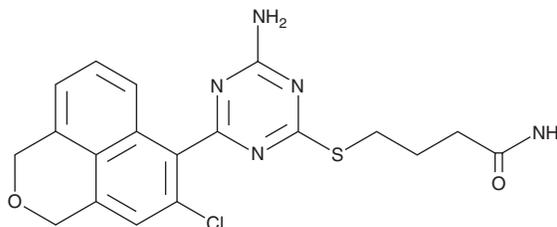
PRODUCT INFORMATION



CH5138303

Item No. 22417

CAS Registry No.: 959763-06-5
Formal Name: 4-[[4-amino-6-(5-chloro-1H,3H-naphtho[1,8-cd]pyran-6-yl)-1,3,5-triazin-2-yl]thio]-butanamide
MF: C₁₉H₁₈ClN₅O₂S
FW: 415.9
Purity: ≥99%
UV/Vis.: λ_{max}: 229 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

CH5138303 is supplied as a crystalline solid. A stock solution may be made by dissolving the CH5138303 in the solvent of choice, which should be purged with an inert gas. CH5138303 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CH5138303 in these solvents is approximately 33 mg/ml.

CH5138303 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CH5138303 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CH5138303 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CH5138303 is a triazine Hsp90 inhibitor.¹ CH5138303 binds to the N-terminus of Hsp90α (K_d = 0.52 nM) and inhibits growth of HCT116 and NCI-N87 human tumor cell lines *in vitro* (IC₅₀s = 89 and 66 nM, respectively). CH5138303 also inhibits growth of human NCI-N87 gastric cancer xenografts (ED₅₀ = 3.9 mg/kg) and displays high oral bioavailability in mice.

Reference

1. Suda, A., Kawasaki, K., Komiyama, S., *et al.* Design and synthesis of 2-amino-6-(1H,3H-benzo[de]isochromen-6-yl)-1,3,5-triazines as novel Hsp90 inhibitors. *Bioorg. Med. Chem.* **22(2)**, 892-905 (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, 10/26/2022

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