

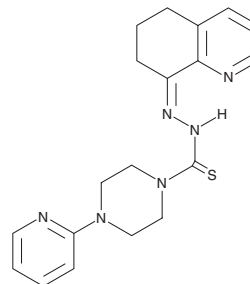
PRODUCT INFORMATION



COTI-2

Item No. 20031

CAS Registry No.: 1039455-84-9
Formal Name: 4-(2-pyridinyl)-2-(6,7-dihydro-8(5H)-quinolinylidene)hydrazide-1-piperazinecarbothioic acid
MF: C₁₉H₂₂N₆S
FW: 366.5
Purity: ≥95% (E/Z mixture)
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

COTI-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the COTI-2 in the solvent of choice. COTI-2 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of COTI-2 in these solvents is approximately 1 and 2.5 mg/ml, respectively.

COTI-2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, COTI-2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. COTI-2 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

COTI-2 is an orally available thiosemicarbazone that activates mutant forms of p53. It induces apoptosis in a wide variety of human tumor cells in culture, inhibits the proliferation of colorectal cancer cell lines more effectively than cetuximab and erlotinib (Item No. 10483), and is active against human glioblastoma cell lines at nanomolar concentrations.¹ COTI-2 functions well *in vivo*, inhibiting the growth of HT-29 colorectal, SHP-77 small cell lung cancer, and OVCAR-3 ovarian carcinoma xenograft cells in mice with minimal toxicity.¹

Reference

1. Salim, K. Y., Vareki, S. M., Danter, W. R., *et al.* COTI-2, a novel small molecule that is active against multiple human cancer cell lines *in vitro* and *in vivo*. *Oncotarget* **7**(27), (2016).

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, 06/28/2022

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