

PRODUCT INFORMATION



CID-5951923

Item No. 21649

CAS Registry No.: 749872-43-3
Formal Name: 3-(3-nitrophenyl)-2-propenoic acid,
2-[methyl(tetrahydro-1,1-dioxido-3-
thienyl)amino]-2-oxoethyl ester

MF: C₁₆H₁₈N₂O₇S

FW: 382.4

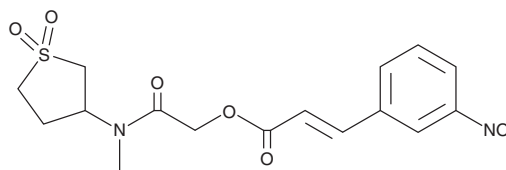
Purity: ≥98%

UV/Vis.: λ_{max}: 260 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

CID-5951923 is supplied as a crystalline solid. A stock solution may be made by dissolving the CID-5951923 in the solvent of choice, which should be purged with an inert gas. CID-5951923 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CID-5951923 in these solvents is approximately 16 and 20 mg/ml, respectively.

CID-5951923 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CID-5951923 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CID-5951923 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CID-5951923 is an inhibitor of the transcription factor Krüppel-like factor 5 (KLF5; IC₅₀ = 2.3 μM).¹ It decreases KLF5 expression in and reduces growth of DLD-1 colorectal adenocarcinoma cells in a concentration-dependent manner. CID-5951923 selectively inhibits proliferation of cancer cell lines that highly express KLF5, particularly colon adenocarcinomas, in the NCI-60 panel of cancer cell lines.

Reference

1. Bialkowska, A.B., Crisp, M., Bannister, T., *et al.* Identification of small-molecule inhibitors of the colorectal cancer oncogene Krüppel-like factor 5 expression by ultrahigh-throughput screening. *Mol. Cancer Ther.* **10(11)**, 2043-2051 (2011).

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, 11/02/2022

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