

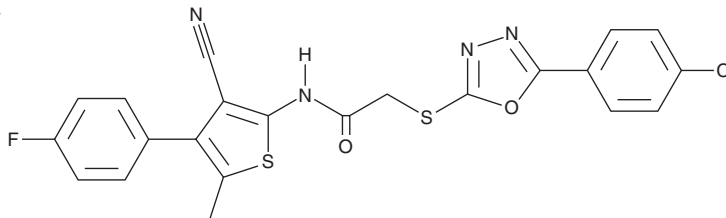
PRODUCT INFORMATION



TG2-179-1

Item No. 38874

CAS Registry No.: 1135023-19-6
Formal Name: 2-[[5-(4-chlorophenyl)-1,3,4-oxadiazol-2-yl]thio]-N-[3-cyano-4-(4-fluorophenyl)-5-methyl-2-thienyl]-acetamide
MF: $C_{22}H_{14}ClFN_4O_2S_2$
FW: 485.0
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 244, 279 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 3 years



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

Laboratory Procedures

TG2-179-1 is supplied as a solid. A stock solution may be made by dissolving the TG2-179-1 in the solvent of choice, which should be purged with an inert gas. TG2-179-1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TG2-179-1 in DMSO is approximately 1 mg/ml and approximately 2 mg/ml in ethanol and DMF.

Description

TG2-179-1 is an inhibitor of BRCA1-associated protein 1 (BAP1).¹ It inhibits the deubiquitinase activity of BAP1 in a cell-free assay when used at concentrations ranging from 25 μM to 1.5 mM. TG2-179-1 is cytotoxic against a panel of eight colon cancer cell lines ($IC_{50}s = 4.48-7.52 \mu M$). *In vivo*, TG2-179-1 (10 and 30 mg/kg) reduces tumor volume in an HCT116 colon cancer mouse xenograft model.

Reference

1. Kang, M., Park, S.G., Lee, S.-A., *et al.* Targeting BAP1 with small compound inhibitor for colon cancer treatment. *Sci. Rep.* **13**(1), 2264 (2023).

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, 07/14/2023

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