

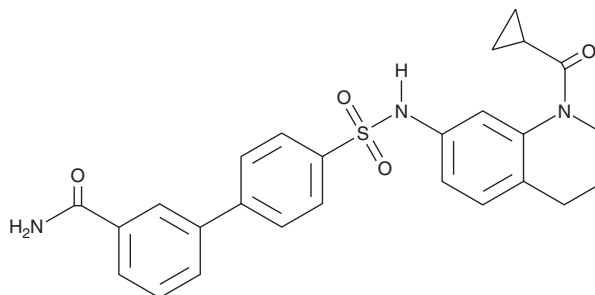
PRODUCT INFORMATION



SU 0268

Item No. 36064

CAS Registry No.: 2210228-45-6
Formal Name: 4'-[[[1-(cyclopropylcarbonyl)-1,2,3,4-tetrahydro-7-quinoliny]amino]sulfonyl]-[1,1'-biphenyl]-3-carboxamide
MF: C₂₆H₂₅N₃O₄S
FW: 475.6
Purity: ≥98%
UV/Vis.: λ_{max}: 254 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

SU 0268 is supplied as a crystalline solid. A stock solution may be made by dissolving the SU 0268 in the solvent of choice, which should be purged with an inert gas. SU 0268 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SU 0268 in ethanol is approximately 15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

SU 0268 is an inhibitor of 8-oxoguanine DNA glycosylase 1 (OGG1; IC₅₀ = 0.059 μM).¹ It is selective for OGG1 over the DNA repair enzymes MTH1, dUTPase, NUDT16, ABH2, ABH3, and SMUG1 at 20 μM. SU 0268 (0.5 μM) inhibits OGG1 and increases the number of 8-oxoguanine DNA lesions in HeLa cell lysates.

Reference

1. Tahara, Y.-K., Auld, D., Ji, D., *et al.* Potent and selective inhibitors of 8-oxoguanine DNA glycosylase. *J. Am. Chem. Soc.* **140**(6), 2105-2114 (2018).

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, 12/19/2022

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