

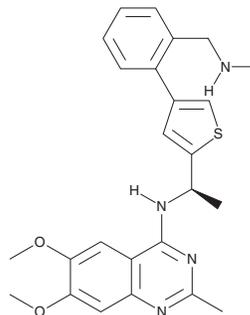
PRODUCT INFORMATION



BAY-293

Item No. 33606

CAS Registry No.: 2244904-70-7
Formal Name: 6,7-dimethoxy-2-methyl-N-[(1R)-1-[4-[2-[(methylamino)methyl]phenyl]-2-thienyl]ethyl]-4-quinazolinamine
MF: C₂₅H₂₈N₄O₂S
FW: 448.6
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 250, 328, 342 nm
Supplied as: A 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

BAY-293 is supplied as a solid. A stock solution may be made by dissolving the BAY-293 in the solvent of choice, which should be purged with an inert gas. BAY-293 is soluble in DMSO.

Description

BAY-293 is an inhibitor of the protein-protein interaction between K-Ras and the guanine nucleotide exchange factor son of sevenless homolog 1 (SOS1; IC₅₀ = 0.021 μM).¹ It inhibits proliferation of K562 and MOLM-13 cells expressing wild-type K-Ras, as well as NCI H358 and Calu-1 cells expressing mutant K-Ras^{G12C} with IC₅₀ values of 1.09, 0.995, 3.48, and 3.19 μM, respectively.

Reference

- Hillig, R.C., Sautier, B., Schroeder, J., *et al.* Discovery of potent SOS1 inhibitors that block RAS activation via disruption of the RAS-SOS1 interaction. *Proc. Natl. Acad. Sci. USA* **116**(7), 2551-2560 (2019).

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, 12/08/2022

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