

# PRODUCT INFORMATION



**BI-2493**

Item No. 44474

**CAS Registry No.:** 2937344-16-4  
**Formal Name:** (4'S)-2'-amino-5,6,6',7'-tetrahydro-3-[2-[(2S)-2-methyl-1-piperazinyl]-4-pyrimidinyl]-spiro[1,2-benzisoxazole-7(4H),4'(5'H)-benzo[b]thiophene]-3'-carbonitrile

**MF:** C<sub>24</sub>H<sub>27</sub>N<sub>7</sub>OS

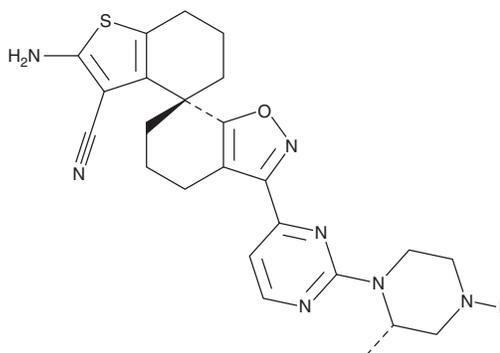
**FW:** 461.6

**Purity:** ≥98%

**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

BI-2493 is supplied as a solid. A stock solution may be made by dissolving the BI-2493 in the solvent of choice, which should be purged with an inert gas. BI-2493 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

## Description

BI-2493 is an inhibitor of mutant K-Ras.<sup>1</sup> It selectively decreases the proliferation of Ba/F3 cells expressing K-Ras containing a glycine-to-cysteine substitution at position 12 (K-Ras<sup>G12C</sup>), K-Ras<sup>G12D</sup>, or K-Ras<sup>G12V</sup> (EC<sub>50</sub>s = 118, 52, and 81 nM, respectively) over Ba/F3 cells expressing H-Ras or N-Ras (EC<sub>50</sub>s = 2,359 and 1,576 nM, respectively). BI-2493 (30 mg/kg twice per day) reduces tumor volume in SW480 colon adenocarcinoma and NCI H358 non-small cell lung cancer (NSCLC) mouse xenograft models.

## Reference

1. Bröker, J., Waterson, A.G., Hodges, T.R., *et al.* Discovery of BI-2493, a pan-KRAS inhibitor showing *in vivo* efficacy. *J. Med. Chem.* **68**(15), 15649-15668 (2025).

### 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, 11/10/2025

## 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