

# PRODUCT INFORMATION



## S49076 (hydrochloride)

Item No. 41173

**CAS Registry No.:** 1265965-19-2  
**Formal Name:** 3-[[[(3Z)-2,3-dihydro-3-[[4-(4-morpholinylmethyl)-1H-pyrrol-2-yl]methylene]-2-oxo-1H-indol-5-yl]methyl]-2,4-thiazolidinedione, monohydrochloride

**MF:** C<sub>22</sub>H<sub>22</sub>N<sub>4</sub>O<sub>4</sub>S • HCl

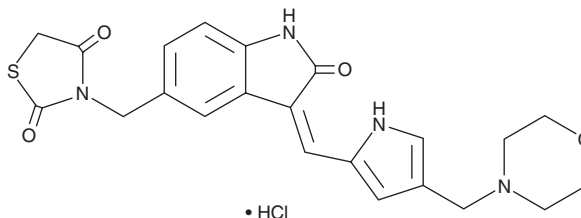
**FW:** 475.0

**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

S49076 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the S49076 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. S49076 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

S49076 is a multi-kinase inhibitor.<sup>1</sup> It inhibits c-Met, Axl, Mer, FGFR1, FGFR2, and FGFR3 (IC<sub>50</sub>s = 1, 7, 2, 18, 17, and 15 nM, respectively). S49076 inhibits Axl- and c-Met-mediated Akt phosphorylation in MDA-MB-231 cells (IC<sub>50</sub>s = 33 and <3 nM, respectively). It also inhibits FGFR1, FGFR2, and FGFR3 autophosphorylation in NCI 1703, SNU-16, and RT-112 cells, respectively (IC<sub>50</sub>s = 50-200 nM). *In vivo*, S49076 (6.25-50 mg/kg) reduces mean tumor volume in GTL-16 and U87MG mouse xenograft models.

### Reference

1. Burbridge, M.F., Bossard, C.J., Saunier, C., *et al.* S49076 is a novel kinase inhibitor of MET, AXL, and FGFR with strong preclinical activity alone and in association with bevacizumab. *Mol. Cancer Ther.* **12(9)**, 1749-1762 (2013).

#### 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, 07/10/2024

#### 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