

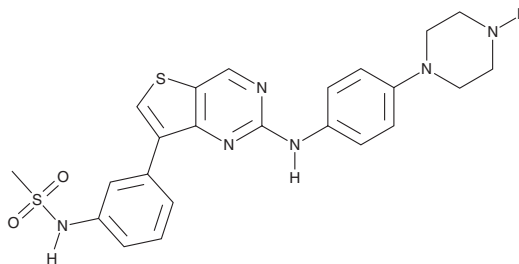
# PRODUCT INFORMATION



## FLT3-IN-17

Item No. 45396

**CAS Registry No.:** 2758999-62-9  
**Formal Name:** N-[3-[2-[[4-(1-piperazinyl)phenyl]amino]thieno[3,2-d]pyrimidin-7-yl]phenyl]-methanesulfonamide  
**MF:** C<sub>23</sub>H<sub>24</sub>N<sub>6</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 480.6  
**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

FLT3-IN-17 is supplied as a solid. A stock solution may be made by dissolving the FLT3-IN-17 in the solvent of choice, which should be purged with an inert gas. FLT3-IN-17 is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

FLT3-IN-17 is a dual inhibitor of FMS-related tyrosine kinase 3 (FLT3) expressing a D835Y mutation associated with quizartinib resistance and focal adhesion kinase (FAK; IC<sub>50</sub>s = <0.5 and 12 nM in cell-free assays).<sup>1</sup> It inhibits the growth of HCT116 colorectal cancer cells, MDA-MB-231 breast cancer cells, and A375 human melanoma cells (GI<sub>50</sub>s = 0.25, 0.46, and 0.49 μM, respectively).

### Reference

1. Cho, H., Shin, I., Yoon, H., *et al.* Identification of thieno[3,2-d]pyrimidine derivatives as dual inhibitors of focal adhesion kinase and FMS-like tyrosine kinase 3. *J. Med. Chem.* **64**(16), 11934-11957 (2021).

#### 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, 04/03/2026

#### 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