

# PRODUCT INFORMATION

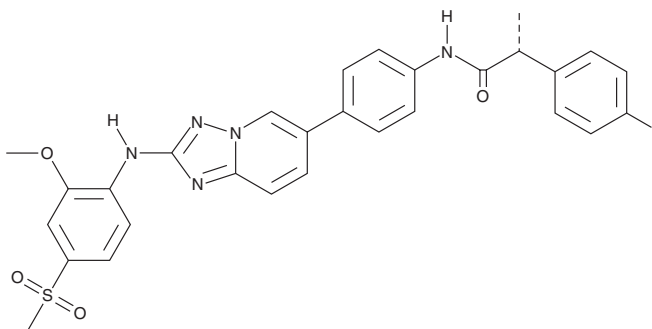


## BAY-1161909

Item No. 41108

**CAS Registry No.:** 1443763-60-7  
**Formal Name:** (αR)-4-fluoro-N-[4-[2-[[2-methoxy-4-(methylsulfonyl)phenyl]amino][1,2,4]triazolo[1,5-a]pyridin-6-yl]phenyl]-α-methyl-benzeneacetamide

**Synonym:** Empesertib  
**MF:** C<sub>29</sub>H<sub>26</sub>FN<sub>5</sub>O<sub>4</sub>S  
**FW:** 559.6  
**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

BAY-1161909 is supplied as a solid. A stock solution may be made by dissolving the BAY-1161909 in the solvent of choice, which should be purged with an inert gas. BAY-1161909 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

### Description

BAY-1161909 is an inhibitor of monopolar spindle 1 (Mps1; IC<sub>50</sub> = 0.34 nM).<sup>1</sup> It inhibits metaphase arrest induced by the antimetabolic agent nocodazole (Item No. 13857) in HeLa cells (IC<sub>50</sub> = 56 nM). BAY-1161909 inhibits proliferation in a panel of 86 cancer cell lines (median IC<sub>50</sub> = 6.7 nM). *In vivo*, BAY-1161909 (2.5 mg/kg) reduces tumor area in a cisplatin-resistant A2780/CP70 ovarian cancer mouse xenograft model.

### Reference

1. Wengner, A.M., Siemeister, G., Koppitz, M., *et al.* Novel Mps1 kinase inhibitors with potent antitumor activity. *Mol. Cancer Ther.* **15**(4), 583-592 (2016).

#### 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, 06/24/2024

#### 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