

# PRODUCT INFORMATION

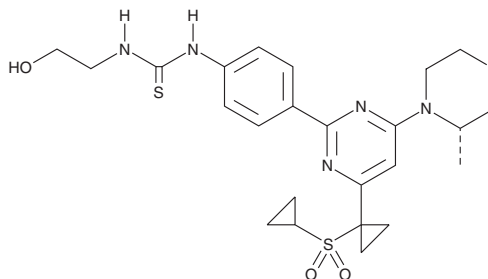


## AZD 3147

Item No. 22474

**CAS Registry No.:** 1101810-02-9  
**Formal Name:** N-[4-[4-[1-(cyclopropylsulfonyl)cyclopropyl]-6-[(3S)-3-methyl-4-morpholinyl]-2-pyrimidinyl]phenyl]-N'-(2-hydroxyethyl)-thiourea

**MF:** C<sub>24</sub>H<sub>31</sub>N<sub>5</sub>O<sub>4</sub>S<sub>2</sub>  
**FW:** 517.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 266, 310 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

AZD 3147 is supplied as a solid. A stock solution may be made by dissolving the AZD 3147 in the solvent of choice, which should be purged with an inert gas. AZD 3147 is soluble in organic solvents such as ethanol, DMSO, and hydrochloric acid. The solubility of AZD 3147 in DMSO and hydrochloric acid is approximately 100 mM and approximately 1 mg/ml in ethanol.

### Description

AZD 3147 is an orally bioavailable dual inhibitor of mammalian target of rapamycin complex 1 (mTORC1) and mTORC2 (IC<sub>50</sub>s = 40.7 and 5.75 nM, respectively, in MDA-MB-468 cells).<sup>1</sup> It is selective for mTOR over PI3Kα, PI3Kβ, PI3Kδ, and PI3Kγ (IC<sub>50</sub>s = 1.51, 912, 5,495, 9,333, and 6,310 nM, respectively, in enzyme assays).

### Reference

1. Pike, K.G., Morris, J., Ruston, L., et al. Discovery of AZD3147: A potent, selective dual inhibitor of mTORC1 and mTORC2. *J. Med. Chem.* **58(5)**, 2326-2349 (2015).

#### 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, 12/15/2022

#### 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