

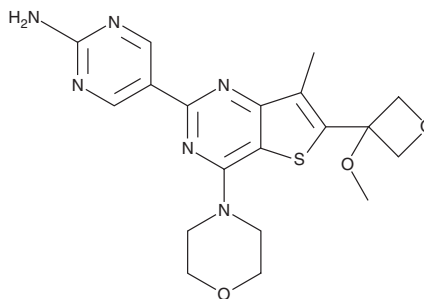
# PRODUCT INFORMATION



## GNE-317

Item No. 27671

**CAS Registry No.:** 1394076-92-6  
**Formal Name:** 5-[6-(3-methoxy-3-oxetanyl)-7-methyl-4-(4-morpholinyl)thieno[3,2-d]pyrimidin-2-yl]-2-pyrimidinamine  
**MF:** C<sub>19</sub>H<sub>22</sub>N<sub>6</sub>O<sub>3</sub>S  
**FW:** 414.5  
**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

GNE-317 is supplied as a solid. A stock solution may be made by dissolving the GNE-317 in the solvent of choice, which should be purged with an inert gas. GNE-317 is soluble in DMSO.

### Description

GNE-317 is a dual inhibitor of PI3K $\alpha$  and the mammalian target of rapamycin (mTOR; apparent K<sub>i</sub>s = 2 and 9 nM, respectively).<sup>1</sup> It inhibits the proliferation of PC3 prostate cancer and A172 glioblastoma cells with EC<sub>50</sub> values of 132 and 240 nM, respectively. GNE-317 (10-40 mg/kg) reduces tumor growth in a U87MG/M glioblastoma mouse xenograft model.

### Reference

1. Heffron, T.P., Salphati, L., Alicke, B., *et al.* The design and identification of brain penetrant inhibitors of phosphoinositide 3-kinase  $\alpha$ . *J. Med. Chem.* **55**(18), 8007-8020 (2012).

#### 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, 11/08/2022

#### 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](http://WWW.CAYMANCHEM.COM)