

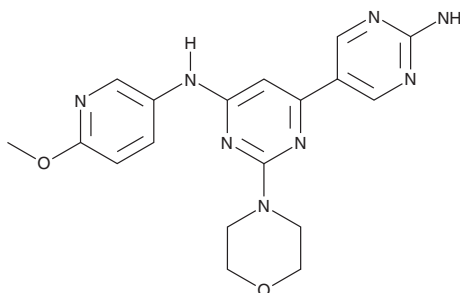
# PRODUCT INFORMATION



## NIBR17

Item No. 37312

**CAS Registry No.:** 944396-88-7  
**Formal Name:** N<sup>6</sup>-(6-methoxy-3-pyridinyl)-2-(4-morpholinyl)-[4,5'-bipyrimidine]-2',6-diamine  
**MF:** C<sub>18</sub>H<sub>20</sub>N<sub>8</sub>O<sub>2</sub>  
**FW:** 380.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 273, 334 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

NIBR17 is supplied as a solid. A stock solution may be made by dissolving the NIBR17 in the solvent of choice, which should be purged with an inert gas. NIBR17 is soluble in the organic solvent DMSO at a concentration of approximately 3 mg/ml.

### Description

NIBR17 is an inhibitor of PI3Kα (IC<sub>50</sub> = <0.002 μM).<sup>1</sup> It inhibits phosphorylation of Akt in, and proliferation of, A2780 ovarian cancer cells (EC<sub>50</sub>s = 0.09 and 0.73 μM, respectively). NIBR17 (60 mg/kg) reduces tumor volume in an A2780 mouse xenograft model.

### Reference

1. Burger, M.T., Knapp, M., Wagman, A.S., *et al.* Synthesis and in vitro and in vivo evaluation of phosphoinositide-3-kinase inhibitors. *ACS Med. Chem. Lett.* **2**(1), 34-38 (2010).

#### 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, 02/08/2023

#### 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)