

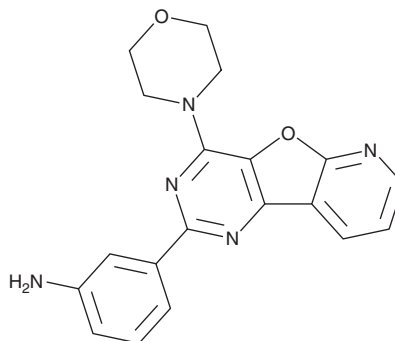
# PRODUCT INFORMATION



**MDK34597**

Item No. 33137

**CAS Registry No.:** 371934-59-7  
**Formal Name:** 3-[4-(4-morpholinyl)pyrido[3',2':4,5]furo[3,2-d]pyrimidin-2-yl]-benzenamine  
**MF:** C<sub>19</sub>H<sub>17</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 347.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232, 285 nm  
**Supplied as:** A crystalline 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

MDK34597 is supplied as a crystalline solid. A stock solution may be made by dissolving the MDK34597 in the solvent of choice, which should be purged with an inert gas. MDK34597 is soluble in DMSO.

## Description

MDK34697 is an inhibitor of PI3K p110α (IC<sub>50</sub> = 141.25 nM).<sup>1</sup>

## Reference

1. Li, Y., Wang, Y., and Zhang, F. Pharmacophore modeling and 3D-QSAR analysis of phosphoinositide 3-kinase p110α inhibitors. *J. Mol. Model* **16(9)**, 1449-1460 (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, 11/14/2022

## 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