

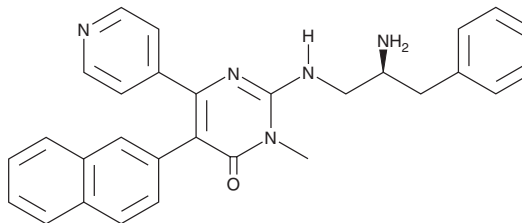
# PRODUCT INFORMATION



## AMG 548

Item No. 34577

**CAS Registry No.:** 864249-60-5  
**Formal Name:** 2-[[[(2S)-2-amino-3-phenylpropyl] amino]-3-methyl-5-(2-naphthalenyl)-6-(4-pyridinyl)-4(3H)-pyrimidinone  
**MF:** C<sub>29</sub>H<sub>27</sub>N<sub>5</sub>O  
**FW:** 461.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222 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

AMG 548 is supplied as a solid. A stock solution may be made by dissolving the AMG 548 in the solvent of choice, which should be purged with an inert gas. AMG 548 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AMG 548 in these solvents is approximately 30 mg/ml.

AMG 548 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AMG 548 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. AMG 548 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

AMG 548 is an inhibitor of p38α MAPK (K<sub>i</sub> = 0.5 nM).<sup>1</sup> It is selective for p38α over p38γ, p38δ, JNK1, JNK2, and JNK3 (K<sub>i</sub>s = 2,600, 4,100, 11,480, 39, and 61 nM, respectively) and is greater than 1,000-fold selective over a panel of 36 additional kinases but also inhibits p38β (K<sub>i</sub> = 3.6 nM). AMG 548 inhibits LPS-induced production of TNF-α and IL-1β in isolated human whole blood (IC<sub>50</sub>s = 3 and 7 nM, respectively).

### Reference

1. Lee, M.R. and Dominguez, C. MAP kinase p38 inhibitors: Clinical results and an intimate look at their interactions with p38α protein. *Curr. Med. Chem.* **12(25)**, 2979-2994 (2005).

#### 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, 12/05/2022

#### 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