

# PRODUCT INFORMATION

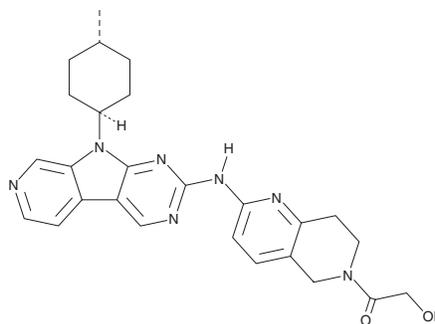


## AMG 925

Item No. 29245

**CAS Registry No.:** 1401033-86-0  
**Formal Name:** 1-[7,8-dihydro-2-[[9-(*trans*-4-methylcyclohexyl)-9H-pyrido[4',3':4,5]pyrrolo[2,3-d]pyrimidin-2-yl]amino]-1,6-naphthyridin-6(5H)-yl]-2-hydroxy-ethanone

**MF:** C<sub>26</sub>H<sub>29</sub>N<sub>7</sub>O<sub>2</sub>  
**FW:** 471.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 227, 268, 342 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

AMG 925 is supplied as a crystalline solid. A stock solution may be made by dissolving the AMG 925 in the solvent of choice, which should be purged with an inert gas. AMG 925 is soluble in the organic solvent chloroform at a concentration of approximately 0.5 mg/ml.

### Description

AMG 925 is a dual inhibitor of FMS-related tyrosine kinase 3 (FLT3) and cyclin-dependent kinase 4 (Cdk4; IC<sub>50</sub>s = 1 and 3 nM, respectively).<sup>1</sup> It is selective for FLT3 and Cdk4 over Cdk1 (IC<sub>50</sub> = 2.22 μM). AMG 925 also binds to FLT3<sup>ITD</sup>, FLT3<sup>D835Y</sup>, FLT3<sup>D835H</sup>, FLT3<sup>K663Q</sup>, and FLT3<sup>N841I</sup> (K<sub>d</sub>s = 1-4 nM). It inhibits proliferation of MOLM-13, COLO 205, and U937 cancer cells (IC<sub>50</sub>s = 19, 55, and 52 nM, respectively). AMG 925 (50, 75, and 150 mg/kg) reduces intratumor levels of pSTAT5 and pRb and inhibits tumor growth in a MOLM-13 acute myeloid leukemia (AML) mouse xenograft model.

### Reference

1. Li, Z., Wang, X., Eksterowicz, J., *et al.* Discovery of AMG 925, a FLT3 and CDK4 dual kinase inhibitor with preferential affinity for the activated state of FLT3. *J. Med. Chem.* **57**(8), 3430-3439 (2014).

#### 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, 10/26/2022

#### 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