

# PRODUCT INFORMATION



## FLT3-IN-3

Item No. 36106

**CAS Registry No.:** 2229050-90-0  
**Formal Name:** N<sup>2</sup>-(*trans*-4-aminocyclohexyl)-9-cyclopentyl-N<sup>6</sup>-[4-(4-morpholinylmethyl)phenyl]-9H-purine-2,6-diamine

**Synonym:** SAN50900

**MF:** C<sub>27</sub>H<sub>38</sub>N<sub>8</sub>O

**FW:** 490.6

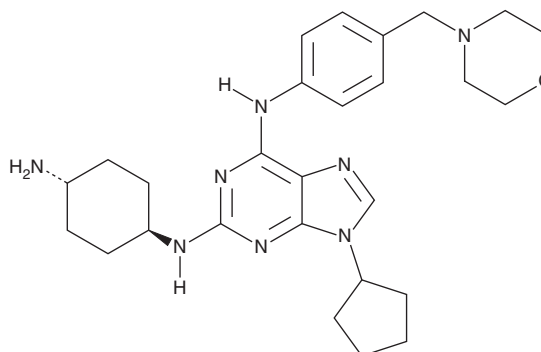
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 262, 312 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

FLT3-IN-3 is supplied as a solid. A stock solution may be made by dissolving the FLT3-IN-3 in the solvent of choice, which should be purged with an inert gas. FLT3-IN-3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of FLT3-IN-3 in ethanol is approximately 15 mg/ml and approximately 10 mg/ml in DMSO and DMF.

### Description

FLT3-IN-3 is an inhibitor of FMS-related tyrosine kinase 3 (FLT3).<sup>1</sup> It inhibits wild-type, FLT3<sup>D835Y</sup>, and FLT3 bearing internal-tandem duplication (FLT3-ITD; IC<sub>50</sub>s = 13, 8, and 3 nM, respectively). FLT3-IN-3 suppresses the growth of MV4-11 acute myeloid leukemia (AML) cells that express FLT3-ITD (GI<sub>50</sub> = 7 nM). It induces cell cycle arrest at the G<sub>1</sub> phase and apoptosis in the same cells. FLT3-IN-3 (10 mg/kg) reduces intratumor FLT3-ITD autophosphorylation in a MV4-11 mouse xenograft model.

### Reference

1. Gucký, T., E., Ř., Radošová Muchová, T., *et al.* Discovery of N<sup>2</sup>-(4-amino-cyclohexyl)-9-cyclopentyl-N<sup>6</sup>-(4-morpholin-4-ylmethyl-phenyl)-9H-purine-2,6-diamine as a potent FLT3 kinase inhibitor for acute myeloid leukemia with FLT3 mutations. *J. Med. Chem.* **61**(90), 3855-3869 (2018).

#### 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, 12/19/2022

#### 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