

# PRODUCT INFORMATION



## ENMD-2076

Item No. 22956

**CAS Registry No.:** 1453868-32-0  
**Formal Name:** 6-(4-methyl-1-piperazinyl)-N-(5-methyl-1H-pyrazol-3-yl)-2-[(1E)-2-phenylethenyl]-4-pyrimidinamine, 2R,3R-dihydroxybutanedioate

**MF:** C<sub>21</sub>H<sub>25</sub>N<sub>7</sub> • C<sub>4</sub>H<sub>6</sub>O<sub>6</sub>  
**FW:** 525.6

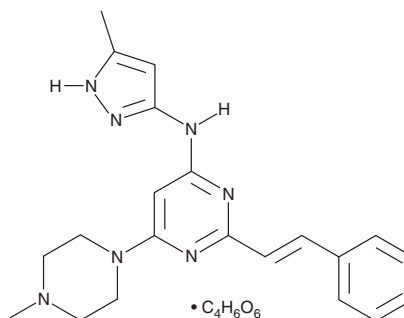
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 203, 269 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

ENMD-2076 is supplied as a crystalline solid. A stock solution may be made by dissolving the ENMD-2076 in the solvent of choice, which should be purged with an inert gas. ENMD-2076 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ENMD-2076 in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

ENMD-2076 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ENMD-2076 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ENMD-2076 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

ENMD-2076 is a multi-kinase inhibitor that inhibits FMS-related tyrosine kinase 3 (FLT3), RET, and Aurora A kinase with IC<sub>50</sub> values of 1.86, 10.4, and 14 nM, respectively.<sup>1</sup> It inhibits additional kinases involved in angiogenesis such as VEGFR3, PDGFRα, and FGFR2, among others, with IC<sub>50</sub> values of less than 100 nM. It inhibits proliferation of several human triple-negative breast cancer (TNBC) cell lines with IC<sub>50</sub> values of less than 1 μM, halts the cell cycle at the G<sub>2</sub> phase, and induces apoptosis.<sup>2</sup> Oral administration of ENMD-2076 (200 mg/kg), reduces proliferation, enhances p53- and p73-mediated cancer cell apoptosis and senescence, and inhibits growth of sensitive primary tumors in a TNBC mouse xenograft model.<sup>3</sup> It also decreases tumor vascular permeability and perfusion and inhibits tumor growth in human TNBC and colorectal cancer mouse xenograft models when administered orally at a dose of 100 mg/kg.<sup>2,4</sup>

### References

1. Fletcher, G.C., Broxk, R.D., Denny, T.A., *et al.* ENMD-2076 is an orally active kinase inhibitor with antiangiogenic and antiproliferative mechanisms of action. *Mol. Cancer Ther.* **10(1)**, 126-137 (2011).
2. Diamond, J.R., Eckhardt, S.G., Tan, A.C., *et al.* Predictive biomarkers of sensitivity to the aurora and angiogenic kinase inhibitor ENMD-2076 in preclinical breast cancer models. *Clin. Cancer Res.* **19(1)**, 291-303 (2013).
3. Ionkina, A.A., Tentler, J.J., Kim, J., *et al.* Efficacy and molecular mechanisms of differentiated response to the aurora and angiogenic kinase inhibitor ENMD-2076 in preclinical models of p53-mutated triple-negative breast cancer. *Front. Oncol.* **7(94)**, (2017).
4. Tentler, J.J., Bradshaw-Pierce, E.L., Serkova, N.J., *et al.* Assessment of the *in vivo* antitumor effects of ENMD-2076, a novel multitargeted kinase inhibitor, against primary and cell line-derived human colorectal cancer xenograft models. *Clin. Cancer Res.* **16(11)**, 2989-2998 (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, 12/02/2022

#### 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