

# PRODUCT INFORMATION



## Golvatinib

Item No. 22143

**CAS Registry No.:** 928037-13-2  
**Formal Name:** N-[2-fluoro-4-[[2-[[[4-(4-methyl-1-piperazinyl)-1-piperidiny]carbonyl]amino]-4-pyridinyl]oxy]phenyl]-N'-(4-fluorophenyl)-1,1-cyclopropanedicarboxamide

**Synonym:** E7050

**MF:** C<sub>33</sub>H<sub>37</sub>F<sub>2</sub>N<sub>7</sub>O<sub>4</sub>

**FW:** 633.7

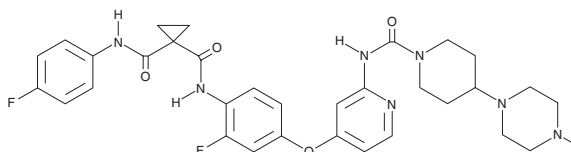
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 244 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

Golvatinib is supplied as a crystalline solid. A stock solution may be made by dissolving the golvatinib in the solvent of choice, which should be purged with an inert gas. Golvatinib is slightly soluble in organic solvents such as DMSO and dimethyl formamide.

### Description

Golvatinib is an orally bioavailable dual inhibitor of the receptor tyrosine kinases c-Met and VEGF receptor 2 (VEGFR2), which are involved in tumor progression.<sup>1</sup> Golvatinib inhibits autophosphorylation of c-Met in MKN45 cells and phosphorylation of VEGFR2 in human umbilical vein endothelial cells (HUVECs; IC<sub>50</sub>s = 14 and 16 nM, respectively). It also inhibits proliferation of a variety of cancer cell lines and of HUVECs stimulated with hepatocyte growth factor (HGF) and VEGF but not bFGF (IC<sub>50</sub>s = 17, 84, and >1,000 nM for HGF-, VEGF-, and bFGF-stimulated HUVECs, respectively). In nude mouse xenograft models using MKN45, Hs746T, SNU-5, or EBC-1 cancer cells, golvatinib (25-200 mg/kg, daily) dose-dependently reduces tumor volume, and it increases survival in the model using MKN45 cells.

### Reference

1. Nakagawa, T., Tohyama, O., Yamaguchi, A., *et al.* E7050: A dual c-Met and VEGFR-2 tyrosine kinase inhibitor promotes tumor regression and prolongs survival in mouse xenograft models. *Cancer Sci.* **101**(1), 210-215 (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](http://WWW.CAYMANCHEM.COM)