

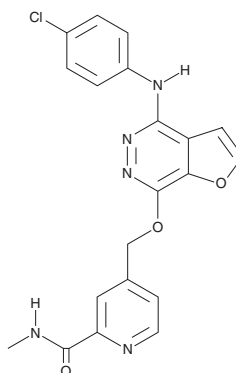
PRODUCT INFORMATION



Telatinib

Item No. 26203

CAS Registry No.: 332012-40-5
Formal Name: 4-[[[4-(4-chlorophenyl)amino]furo[2,3-d]pyridazin-7-yl]oxy)methyl]-N-methyl-2-pyridinecarboxamide
Synonym: BAY 57-9352
MF: C₂₀H₁₆ClN₅O₃
FW: 409.8
Purity: ≥98%
UV/Vis.: λ_{max}: 286 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

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

Telatinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, telatinib should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Telatinib has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Telatinib is a multi-kinase inhibitor that inhibits VEGF receptor 2 (VEGFR2), VEGFR3, PDGFRα, and c-Kit (IC₅₀s = 6, 4, 15, and 1 nM, respectively).¹ It also binds to the transmembrane region of the ABCG2 efflux transporter and enhances intracellular accumulation of [³H]-mitoxantrone in ABCG2-overexpressing cells.² Telatinib (15 mg/kg) decreases tumor growth rate and size in an H460/MX20 mouse xenograft model.

References

1. Steeghs, N., Gelderblom, H., Roodt, J.O., *et al.* Hypertension and rarefaction during treatment with telatinib, a small molecule angiogenesis inhibitor. *Clin. Cancer Res.* **14**(11), 3470-3476 (2008).
2. Sodani, K., Patel, A., Anreddy, N., *et al.* Telatinib reverses chemotherapeutic multidrug resistance mediated by ABCG2 efflux transporter *in vitro* and *in vivo*. *Biochem. Pharmacol.* **89**(1), 52-61 (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, 12/07/2022

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