

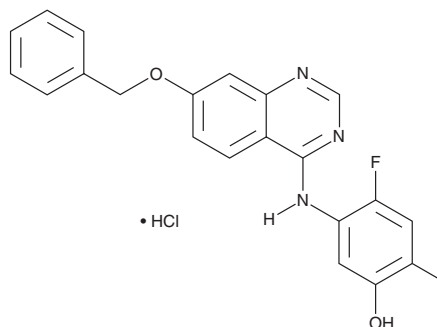
# PRODUCT INFORMATION



## ZM 323881 (hydrochloride)

Item No. 21347

**CAS Registry No.:** 193000-39-4  
**Formal Name:** 4-fluoro-2-methyl-5-[[7-(phenylmethoxy)-4-quinazoliny]amino]-phenol, monohydrochloride  
**MF:** C<sub>22</sub>H<sub>18</sub>FN<sub>3</sub>O<sub>2</sub> • HCl  
**FW:** 411.9  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250, 321 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

ZM 323881 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ZM 323881 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ZM 323881 (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ZM 323881 (hydrochloride) in these solvents is approximately 30 mg/ml.

ZM 323881 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ZM 323881 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ZM 323881 (hydrochloride) 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

ZM 323881 is a potent inhibitor of vascular endothelial growth factor receptor 2 (VEGFR2; IC<sub>50</sub> = <2 nM), a receptor tyrosine kinase.<sup>1</sup> It is selective for VEGFR2 over VEGFR1, PDGFRβ, FGFR1, EGFR, and ErbB2 at concentrations greater than 50 μM. It inhibits VEGF-A-stimulated cell proliferation of human umbilical vein endothelial cells (HUVECs; IC<sub>50</sub> = 8 nM) but not basal proliferation or that stimulated by EGF or bFGF (IC<sub>50</sub>s = >5,000, 1,900, and 1,600 nM, respectively).

### Reference

1. Whittles, C.E., Pocock, T.M., Wedge, S.R., *et al.* ZM323881, a novel inhibitor of vascular endothelial growth factor-receptor-2 tyrosine kinase activity. *Microcirculation* **9**(6), 513-522 (2002).

#### 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