

PRODUCT INFORMATION

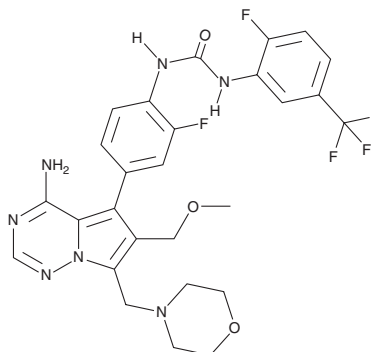


ACTB-1003

Item No. 30069

CAS Registry No.: 939805-30-8
Formal Name: N-[4-[4-amino-6-(methoxymethyl)-7-(4-morpholinylmethyl)pyrrolo[2,1-f][1,2,4]triazin-5-yl]-2-fluorophenyl]-N'-[2-fluoro-5-(trifluoromethyl)phenyl]-urea

MF: C₂₇H₂₆F₅N₇O₃
FW: 591.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ACTB-1003 is supplied as a crystalline solid. A stock solution may be made by dissolving the ACTB-1003 in the solvent of choice, which should be purged with an inert gas. ACTB-1003 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ACTB-1003 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

ACTB-1003 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ACTB-1003 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ACTB-1003 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

ACTB-1003 is a multi-kinase inhibitor (IC₅₀s = 6, 2, 4, 5, and 32 nM for FGFR1, VEGFR2, Tie-2, RSK, and p70S6K, respectively).¹ It inhibits growth of OPM2 human multiple myeloma and murine Ba/F3-TEL-FGFR1 leukemia cells *in vitro*. ACTB-1003 inhibits tumor angiogenesis in an H460 mouse xenograft model. It also reduces tumor growth in an HCT116 mouse xenograft model when administered in combination with 5-fluorouracil (5-FU; Item No. 14416) or paclitaxel (Item No. 10461).

Reference

1. Patel, K., Fattaey, A., and Burd, A. ACTB-1003: An oral kinase inhibitor targeting cancer mutations (FGFR), angiogenesis (VEGFR2, TEK), and induction of apoptosis (RSK and p70S6K). *J. Clin. Oncol.* **28(15_suppl)**, e13665-e13665 (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, 05/11/2020

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