

PRODUCT INFORMATION

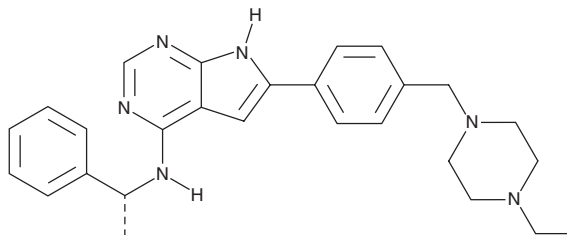


AEE788

Item No. 18416

CAS Registry No.: 497839-62-0
Formal Name: 6-[4-[(4-ethyl-1-piperazinyl)methyl]phenyl]-N-[(1R)-1-phenylethyl]-7H-pyrrolo[2,3-d]pyrimidin-4-amine
Synonym: NVP-AEE788

MF: C₂₇H₃₂N₆
FW: 440.6
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 327 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

AEE788 is supplied as a crystalline solid. A stock solution may be made by dissolving the AEE788 in the solvent of choice, which should be purged with an inert gas. AEE788 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AEE788 in ethanol is approximately 5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

AEE788 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AEE788 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AEE788 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AEE788 is an inhibitor of both epidermal growth factor (EGF) and vascular endothelial growth factor (VEGF) receptor tyrosine kinase family members.¹ It can inhibit EGFR and VEGFR tyrosine kinases with IC₅₀ values of 2, 6, 77, and 59 nM for EGFR, ErbB2, KDR, and FLT1, respectively.¹ AEE788 demonstrates antiproliferative activity against a range of EGFR and ErbB2-overexpressing cell lines and can inhibit the proliferation of EGF- and VEGF-stimulated human umbilical vein endothelial cells.¹ It also has been shown to have antitumor and anti-angiogenic activity in numerous animal models of cancer.¹

Reference

1. Traxler, P., Allegrini, P.R., Brandt, R., et al. AEE788: A dual family epidermal growth factor receptor/ErbB2 and vascular endothelial growth factor receptor tyrosine kinase inhibitor with antitumor and antiangiogenic activity. *Cancer Res.* **64**(14), 4931-4941 (2004).

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, 12/06/2022

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