

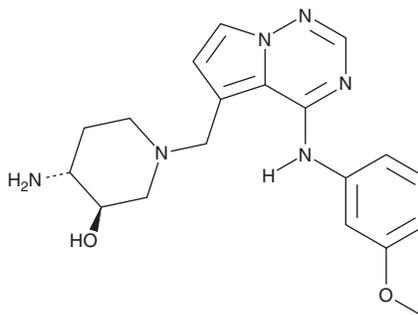
PRODUCT INFORMATION



BMS 690514

Item No. 44552

CAS Registry No.: 859853-30-8
Formal Name: (3R,4R)-4-amino-1-[[4-[(3-methoxyphenyl)amino]pyrrolo[2,1-f][1,2,4]triazin-5-yl)methyl]-3-piperidinol
MF: C₁₉H₂₄N₆O₂
FW: 368.4
Purity: ≥98%
Supplied as: A 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

BMS 690514 is supplied as a solid. A stock solution may be made by dissolving the BMS 690514 in the solvent of choice, which should be purged with an inert gas. BMS 690514 is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

BMS 690514 is a pan-inhibitor of HER and VEGFR family kinases (IC₅₀s = 5, 19, 60, 50, 50, and 25 nM for EGFR, HER2, HER4, VEGFR1, VEGFR2, and VEGFR3, respectively).¹ It is selective for these kinases over 11 additional kinases but also inhibits FMS-related tyrosine kinase 3 (FLT3), calcium/calmodulin-dependent protein kinase II (CAMKII), LCK, and PKA (IC₅₀s = 110, 170, 220, and 280 nM, respectively). BMS 690514 selectively inhibits the proliferation of cancer cells containing EGFR gene amplifications or mutations and HER2 gene amplifications over those expressing wild-type EGFR or not expressing EGFR and HER2. It inhibits tumor growth and angiogenesis in an L2987 lung cancer mouse xenograft model when administered at a dose of 90 mg/kg.

Reference

1. Wong, T.W., Lee, F.Y., Emanuel, S., *et al.* Antitumor and antiangiogenic activities of BMS-690514, an inhibitor of human EGF and VEGF receptor kinase families. *Clin. Cancer Res.* **17(12)**, 4031-4041 (2011).

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, 12/19/2025

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