

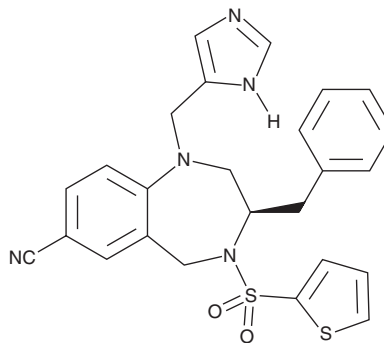
PRODUCT INFORMATION



BMS 214662

Item No. 28713

CAS Registry No.: 195987-41-8
Formal Name: (3R)-2,3,4,5-tetrahydro-1-(1H-imidazol-5-ylmethyl)-3-(phenylmethyl)-4-(2-thienylsulfonyl)-1H-1,4-benzodiazepine-7-carbonitrile
MF: C₂₅H₂₃N₅O₂S₂
FW: 489.6
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 214662 is supplied as a solid. A stock solution may be made by dissolving the BMS 214662 in the solvent of choice, which should be purged with an inert gas. BMS 214662 is slightly soluble in methanol and DMSO.

Description

BMS 214662 is a potent inhibitor of farnesyltransferase (FTase; IC₅₀ = 1.3 nM).¹ It is selective for FTase over geranylgeranyl transferase (GGTase; IC₅₀ = 1,900 nM). It inhibits the growth of MEK2, A2780, and PC3 cancer cells expressing wild-type Ras (IC₅₀s = 2.5, 0.04, and 0.15 μM, respectively), as well as HCT116, MIP, RC-165, and MIA PaCa-2 cells expressing mutant K-Ras (IC₅₀s = 0.06, 0.3, 0.3, and 0.12 μM, respectively). BMS 214662 induces apoptosis in HCT116 cells in a concentration-dependent manner. *In vivo*, BMS 214662 (600 mg/kg) is curative in an HCT116 mouse xenograft model. It also reduces tumor growth in Calu-1, HT-29, EJ-1, and MIA PaCa-2 mouse xenograft models.

Reference

1. Rose, W.C., Lee, F.Y., Fairchild, C.R., *et al.* Preclinical antitumor activity of BMS-214662, a highly apoptotic and novel farnesyltransferase inhibitor. *Cancer Res.* **61(20)**, 7507-7517 (2001).

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, 11/04/2022

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