

PRODUCT INFORMATION



BGG463

Item No. 41146

CAS Registry No.: 890129-26-7

Formal Name: 6-[[6-(acetylamino)-4-pyrimidinyl]oxy]-N-[4-[(4-methyl-1-piperazinyl)methyl]-3-(trifluoromethyl)phenyl]-1-naphthalenecarboxamide

Synonyms: K03859, NVP-BGG463

MF: $C_{30}H_{29}F_3N_6O_3$

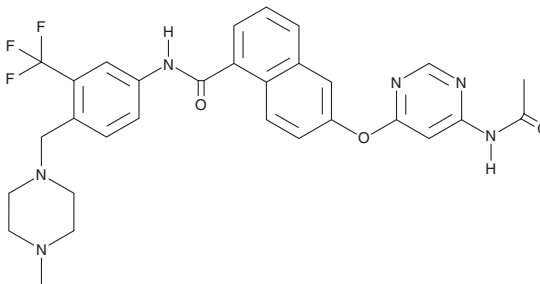
FW: 578.6

Purity: $\geq 95\%$

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

BGG463 is supplied as a solid. A stock solution may be made by dissolving the BGG463 in the solvent of choice, which should be purged with an inert gas. BGG463 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (1-10 mg/ml) in methanol.

Description

BGG463 is a derivative of BBT594, an inhibitor of the fusion protein Bcr-Abi^{T315I}.¹

Reference

1. Manley, P.W., Brügger, J., Floersheimer, A., *et al.* Rational design of T3151 BCR-abl inhibitors for the treatment of drug-resistant chronic myelogenous leukemia (CML): BGG463. *Med. Chem.* **62**(7/8), 579 (2008).

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, 06/07/2024

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