

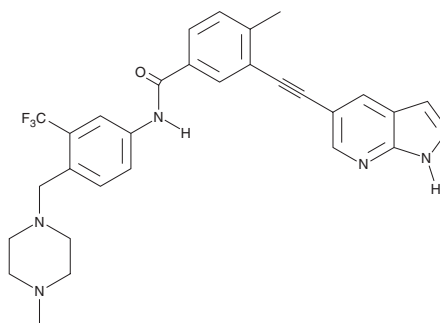
PRODUCT INFORMATION



GZD-824

Item No. 21508

CAS Registry No.: 1257628-77-5
Formal Name: 4-methyl-N-[4-[(4-methyl-1-piperazinyl)methyl]-3-(trifluoromethyl)phenyl]-3-[2-(1H-pyrazolo[3,4-b]pyridin-5-yl)ethynyl]-benzamide
MF: C₂₉H₂₇F₃N₆O
FW: 532.6
Purity: ≥98%
UV/Vis.: λ_{max}: 274, 324 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

GZD-824 is supplied as a crystalline solid. A stock solution may be made by dissolving the GZD-824 in the solvent of choice. GZD-824 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

GZD-824 is an orally available inhibitor of a broad spectrum of Bcr/Abl tyrosine kinase mutants including T315I (IC₅₀s = 0.34 and 0.68 nM for wild-type Bcr/Abl and Bcr/Abl^{T315I}, respectively).¹ It has been shown to suppress the proliferation of Bcr/Abl-positive K562 and Ku812 human chronic myelogenous leukemia cells (IC₅₀s = 0.2 and 0.13 nM, respectively) and induce tumor regression in mouse xenograft tumor models driven by either wild-type or mutant Bcr/Abl.¹

Reference

1. Ren, X., Pan, X., Zhang, Z., *et al.* Identification of GZD824 as an orally bioavailable inhibitor that targets phosphorylated and nonphosphorylated breakpoint cluster region-Abelson (Bcr-Abl) kinase and overcomes clinically acquired mutation-induced resistance against imatinib. *J. Med. Chem.* **56**(3), 879-894 (2013).

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, 10/25/2022

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