

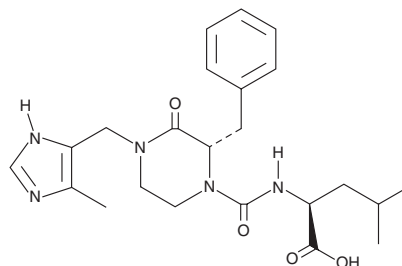
PRODUCT INFORMATION



GGTI 2418

Item No. 28714

CAS Registry No.: 501010-06-6
Formal Name: N-[[[(2S)-4-[(4-methyl-1H-imidazol-5-yl)methyl]-3-oxo-2-(phenylmethyl)-1-piperazinyl]carbonyl]-L-leucine
MF: C₂₃H₃₁N₅O₄
FW: 441.5
Purity: ≥98%
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

GGTI 2418 is supplied as a crystalline solid. A stock solution may be made by dissolving the GGTI 2418 in the solvent of choice, which should be purged with an inert gas. GGTI 2418 is soluble in DMSO.

Description

GGTI 2418 is an inhibitor of geranylgeranyl transferase type I (GGTase I; IC₅₀ = 9.5 nM).¹ It is selective for GGTase I over farnesyltransferase (FTase; IC₅₀ = 53,000 nM). GGTI 2418 (100 mg/kg) reduces tumor growth in an MDA-MB-231 breast cancer mouse xenograft model.

Reference

1. Kazi, A., Carie, A., Blaskovich, M.A., et al. Blockade of protein geranylgeranylation inhibits Cdk2-dependent p27^{Kip1} phosphorylation on Thr187 and accumulates p27^{Kip1} in the nucleus: Implications for breast cancer therapy. *Mol. Cell Biol.* **29**(8), 2254-2263 (2009).

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, 12/08/2022

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