

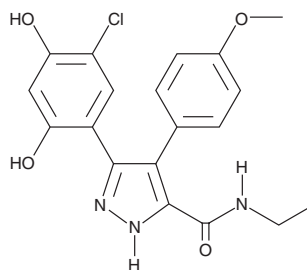
PRODUCT INFORMATION



VER-49009

Item No. 13131

CAS Registry No.: 558640-51-0
Formal Name: 3-(5-chloro-2,4-dihydroxyphenyl)-N-ethyl-4-(4-methoxyphenyl)-1H-pyrazole-5-carboxamide
MF: C₁₉H₁₈ClN₃O₄
FW: 387.8
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.

Description

The heat shock proteins (Hsps) act as molecular chaperones. Hsp90 is an abundant protein with roles in protein folding, cell signaling, and cancer. VER-49009 is a pyrazole compound that inhibits Hsp90 with an IC₅₀ value of 47 nM.¹ It produces a cellular antiproliferative GI₅₀ value of 685 nM when tested against a human cancer cell line panel.¹ Consistent with inhibition of Hsp90, VER-49009 induces the expression of Hsp27 and Hsp72 while reducing the client proteins c-Raf, B-Raf, survivin, and PRMT5, causing cell cycle arrest and apoptosis.¹

Reference

1. Sharp, S.Y., Prodromou, C., Boxall, K., *et al.* Inhibition of the heat shock protein 90 molecular chaperone *in vitro* and *in vivo* by novel, synthetic, potent resorcinylic pyrazole/isoxazole amide analogues. *Mol. Cancer Ther.* **6(4)**, 1198-1211 (2007).

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, 03/11/2024

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