

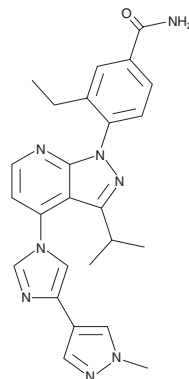
PRODUCT INFORMATION



TAS 116

Item No. 33779

CAS Registry No.: 1260533-36-5
Formal Name: 3-ethyl-4-[3-(1-methylethyl)-4-[4-(1-methyl-1H-pyrazol-4-yl)-1H-imidazol-1-yl]-1H-pyrazolo[3,4-b]pyridin-1-yl]-benzamide
MF: C₂₅H₂₆N₈O
FW: 454.5
Purity: ≥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

TAS 116 is supplied as a solid. A stock solution may be made by dissolving the TAS 116 in the solvent of choice, which should be purged with an inert gas. TAS 116 is soluble in DMSO.

Description

TAS 116 is a heat shock protein 90 (Hsp90) inhibitor.¹ It selectively binds to Hsp90α and Hsp90β (K_is = 34.7 and 21.3 nM, respectively) over glucose-regulated protein 94 (GRP94) and TGF-β receptor-associated protein 1 (TRAP1; K_is = 50,000 nM for both). TAS 116 (3.6, 7.1, and 14 mg/kg) reduces intratumoral protein levels of HER2, HER3, and Akt, as well as reduces tumor growth in a HER2-expressing NCI N87 gastric cancer mouse xenograft model and does not induce photoreceptor injury in rats when administered at a dose of 12 mg/kg per day.

Reference

1. Ohkubo, S., Kodama, Y., Muraoka, H., *et al.* TAS-116, a highly selective inhibitor of heat shock protein 90α and β, demonstrates potent antitumor activity and minimal ocular toxicity in preclinical models. *Mol. Cancer Ther.* **14**(1), 14-22 (2015).

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