

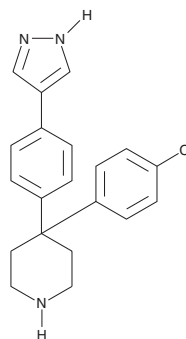
PRODUCT INFORMATION



AT7867

Item No. 16848

CAS Registry No.: 857531-00-1
Formal Name: 4-(4-chlorophenyl)-4-[4-(1H-pyrazol-4-yl)phenyl]-piperidine
MF: C₂₀H₂₀ClN₃
FW: 337.9
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

AT7867 is supplied as a crystalline solid. A stock solution may be made by dissolving the AT7867 in the solvent of choice, which should be purged with an inert gas. AT7867 is soluble in the organic solvent DMSO at a concentration of approximately 0.33 mg/ml.

Description

AT7867 is a potent and orally bioavailable inhibitor of Akt isoforms Akt1, 2, and 3 (IC₅₀s = 32, 17, and 47 nM, respectively).¹ It also inhibits p70S6 kinase and PKA (IC₅₀s = 85 and 20 nM, respectively), but is without effect against a panel of other kinases.¹ AT7867 inhibits growth and induces apoptosis in a variety of cancer cell lines *in vitro* and suppresses tumor growth of PTEN-deficient xenografts in mice.¹

Reference

1. Grimshaw, K.M., Hunter, L.-J.K., Yap, T.A., *et al.* AT7867 is a potent and oral inhibitor of AKT and p70S6 kinase that induces pharmacodynamic changes and inhibits human tumor xenograft growth. *Mol. Cancer Ther.* **9(5)**, 1100-1110 (2010).

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, 10/17/2022

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