

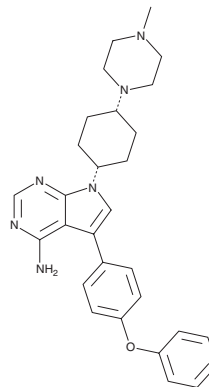
PRODUCT INFORMATION



RK-20448

Item No. 25606

CAS Registry No.: 479501-40-1
Formal Name: 7-[*cis*-4-(4-methyl-1-piperazinyl)cyclohexyl]-5-(4-phenoxyphenyl)-7H-pyrrolo[2,3-*d*]pyrimidin-4-amine
MF: C₂₉H₃₄N₆O
FW: 482.6
Purity: ≥98%
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

RK-20448 is supplied as a solid. A stock solution may be made by dissolving the RK-20448 in the solvent of choice, which should be purged with an inert gas. RK-20448 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

RK-20448 is an ATP-competitive inhibitor of Lck, Src, KDR/VEGF2R, and Tie-2 (IC₅₀s = 0.24, 1.19, 10.74, and 5.85 μM, respectively).¹ It also inhibits BLK, Csk, Fyn, and Lyn (IC₅₀s = 0.37, 4.27, 2.03, and 0.43 μM, respectively). RK-20448 is the *cis* isomer of A-419259 (Item No. 18168).

Reference

1. Calderwood, D.J., Johnston, D.N., Munschauer, R., *et al.* Pyrrolo[2,3-*d*]pyrimidines containing diverse N-7 substituents as potent inhibitors of Lck. *Bioorg. Med. Chem. Lett.* **12(12)**, 1683-1686 (2002).

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, 01/03/2023

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