

# PRODUCT INFORMATION

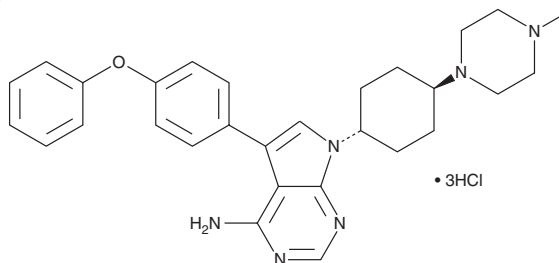


## A-419259 (hydrochloride)

Item No. 18168

**CAS Registry No.:** 1435934-25-0  
**Formal Name:** 7-[*trans*-4-(4-methyl-1-piperazinyl)cyclohexyl]-5-(4-phenoxyphenyl)-7H-pyrrolo[2,3-d]pyrimidin-4-amine, trihydrochloride

**Synonym:** RK-20449  
**MF:** C<sub>29</sub>H<sub>34</sub>N<sub>6</sub>O • 3HCl  
**FW:** 592.0  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 238, 291 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

A-419259 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the A-419259 (hydrochloride) in the solvent of choice. A-419259 (hydrochloride) is soluble in the organic solvent methanol, which should be purged with an inert gas, at a concentration of approximately 1.4 mg/ml.

### Description

A-419259 is an inhibitor of Src family kinases, including Src, LCK, Lyn, and Hck (IC<sub>50</sub>s = 9, <3, <3, and 11.26 nM, respectively).<sup>1,2</sup> It is selective for these kinases over c-Abl (IC<sub>50</sub> = 3,000 nM) and PKC (IC<sub>50</sub> = >33 μM).<sup>1</sup> A-419259 inhibits growth of Philadelphia chromosome-positive (Ph<sup>+</sup>) K-562 and Meg-01 myeloid leukemia cells (IC<sub>50</sub>s = 0.1-0.3 and 0.1 μM, respectively), but not Ph<sup>-</sup> TF-1 and HEL cells. It induces apoptosis in K-562 cells in a concentration-dependent manner. A-419259 (300 nM) inhibits differentiation of murine embryonic stem cells while maintaining pluripotency.<sup>3</sup> It reduces the total number of acute myeloid leukemia (AML) cells, as well as AML stem cells, in the bone marrow and spleen in mouse patient-derived xenograft (PDX) models of AML when administered at a dose of 30 mg/kg twice daily.<sup>4</sup>

### References

1. Wilson, M.B., Schreiner, S.J., Choi, H.J., *et al.* Selective pyrrolo-pyrimidine inhibitors reveal a necessary role for Src family kinases in Bcr-Abl signal transduction and oncogenesis. *Oncogene* **21(53)**, 8075-8088 (2002).
2. Pene-Dumitrescu, T., Peterson, L.F., Donato, N.J., *et al.* An inhibitor-resistant mutant of Hck protects CML cells against the antiproliferative and apoptotic effects of the broad-spectrum Src family kinase inhibitor A-419259. *Oncogene* **27(56)**, 7055-7069 (2008).
3. Meyn, M.A., III, Schreiner, S.J., Dumitrescu, T.P., *et al.* SRC family kinase activity is required for murine embryonic stem cell growth and differentiation. *Mol. Pharmacol.* **68(5)**, 1320-1330 (2005).
4. Saito, Y., Yuki, H., Kuratani, M., *et al.* A pyrrolo-pyrimidine derivative targets human primary AML stem cells *in vivo*. *Sci. Transl. Med.* **5(181)**:181ra52, (2013).

#### 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, 08/02/2018

#### 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