

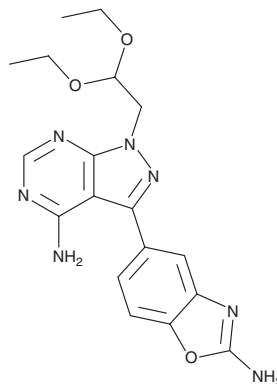
PRODUCT INFORMATION



eCF309

Item No. 21118

CAS Registry No.: 2001571-40-8
Formal Name: 3-(2-amino-5-benzoxazolyl)-
1-(2,2-diethoxyethyl)-1H-
pyrazolo[3,4-d]pyrimidin-4-amine
MF: C₁₈H₂₁N₇O₃
FW: 383.4
Purity: ≥98%
UV/Vis.: λ_{max}: 242, 288, 411 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

eCF309 is supplied as a crystalline solid. A stock solution may be made by dissolving the eCF309 in the solvent of choice, which should be purged with an inert gas. eCF309 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of eCF309 in these solvents is approximately 2 and 1.6 mg/ml, respectively.

eCF309 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, eCF309 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. eCF309 has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

eCF309 is an ATP-competitive inhibitor of mTOR (IC₅₀ = 15 nM *in vitro* and in cells) with >65-fold selectivity over other kinases, including PI3Ks.¹

Reference

1. Fraser, C., Carragher, N.O., and Unciti-Broceta, A. eCF309: A potent, selective and cell-permeable mTOR inhibitor. *Med. Chem. Commun.* **7**(3), 471-477 (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, 10/07/2022

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