

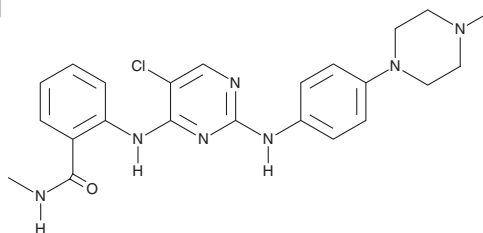
PRODUCT INFORMATION



CTX-0294885

Item No. 22197

CAS Registry No.: 1439934-41-4
Formal Name: 2-[[5-chloro-2-[[4-(1-piperazinyl)phenyl]amino]-4-pyrimidinyl]amino]-N-methylbenzamide
MF: C₂₂H₂₄ClN₇O
FW: 437.9
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 286 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

CTX-0294885 is supplied as a crystalline solid. A stock solution may be made by dissolving the CTX-0294885 in the solvent of choice, which should be purged with an inert gas. CTX-0294885 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CTX-0294885 in ethanol is approximately 0.1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

CTX-0294885 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CTX-0294885 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CTX-0294885 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CTX-0294885 is a broad spectrum kinase inhibitor that inhibited FAK, FLT3, JAK2, JAK3, Src, Aurora kinase A, and VEGF receptor 3 (IC₅₀s = 4, 1, 3, 28, 2, 18, and 3 nM, respectively) in an initial screen.¹ In an affinity purification experiment, CTX-0294885 beads bound to kinases from every major group and captured all members of the AKT family of proteins. In a large-scale kinome profiling experiment, it captured 235 kinases from MDA-MB-231 breast cancer cells.

Reference

1. Zhang, L., Holmes, I.P., Hochgräfe, F., *et al.* Characterization of the novel broad-spectrum kinase inhibitor CTX-0294885 as an affinity reagent for mass spectrometry-based kinome profiling. *J. Proteome Res.* **12**(7), 3104-3116 (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, 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