

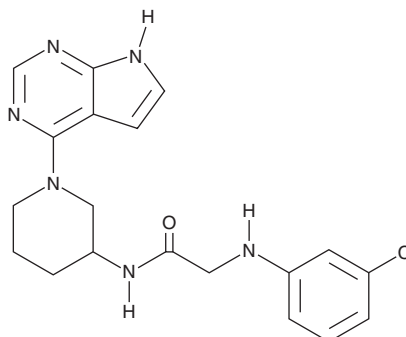
PRODUCT INFORMATION



SNS-062 analog

Item No. 29200

CAS Registry No.: 1270014-40-8
Formal Name: 2-[(3-chlorophenyl)amino]-N-[1-(7H-pyrrolo[2,3-d]pyrimidin-4-yl)-3-piperidinyl]-acetamide
Synonym: BTK-IN-1
MF: C₁₉H₂₁ClN₆O
FW: 384.9
Purity: ≥98%
UV/Vis.: λ_{max}: 252, 278 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

SNS-062 analog is supplied as a crystalline solid. A stock solution may be made by dissolving the SNS-062 analog in the solvent of choice, which should be purged with an inert gas. SNS-062 analog is soluble in the organic solvent DMSO.

Description

SNS-062 analog is an inhibitor of Bruton's tyrosine kinase (BTK; IC₅₀ = <100 nM).¹

Reference

1. Bui, M., Conlon, P., Erlanson, D.A., *et al.* Heteroaryl BTK inhibitors. **WO 2011/029043 A1** (2011).

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, 10/21/2022

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