

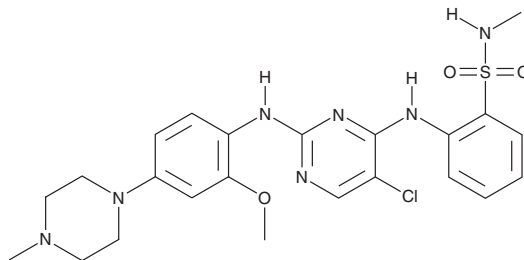
PRODUCT INFORMATION

ALK Inhibitor 2

Item No. 36823

CAS Registry No.: 761438-38-4
Formal Name: 2-[[5-chloro-2-[[2-methoxy-4-(4-methyl-1-piperazinyl)phenyl]amino]-4-pyrimidinyl]amino]-N-methyl-benzenesulfonamide

MF: C₂₃H₂₈ClN₇O₃S
FW: 518.0
Purity: ≥98%
UV/Vis.: λ_{max}: 287 nm
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

ALK inhibitor 2 is supplied as a solid. A stock solution may be made by dissolving the ALK inhibitor 2 in the solvent of choice, which should be purged with an inert gas. ALK inhibitor 2 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ALK inhibitor 2 in these solvents is approximately 2 and 1 mg/ml, respectively.

Description

ALK inhibitor 2 is an inhibitor of testis-specific serine kinase 2 (TSSK2; IC₅₀ = 37 nM).¹

Reference

1. Hawkinson, J.E., Sinville, R., Mudaliar, D., *et al.* Potent pyrimidine and pyrrolopyrimidine inhibitors of testis-specific serine/threonine kinase 2 (TSSK2). *ChemMedChem* **12**(22), 1857-1865 (2017).

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, 12/21/2022

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