

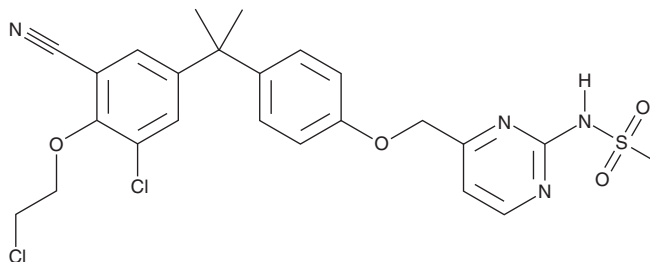
PRODUCT INFORMATION



Masofaniten Item No. 41670

CAS Registry No.: 2416716-62-4
Formal Name: N-[4-[[4-[1-[3-chloro-4-(2-chloroethoxy)-5-cyanophenyl]-1-methylethyl]phenoxy]methyl]-2-pyrimidinyl]-methanesulfonamide

Synonym: EPI-7386
MF: C₂₄H₂₄Cl₂N₄O₄S
FW: 535.4
Purity: ≥95%
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

Masofaniten is supplied as a solid. A stock solution may be made by dissolving the masofaniten in the solvent of choice, which should be purged with an inert gas. Masofaniten is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Masofaniten is an androgen receptor N-terminal transactivation domain inhibitor (ARTADI).¹ It binds to the tryptophan residues in positions 397 (W³⁹⁷) and W⁴³³ in the AR TAD.

Reference

1. Wang, C., Hockerman, S., Jacobsen, E.J., *et al.* Selective inhibition of the p38α MAPK-MK2 axis inhibits inflammatory cues including inflammasome priming signals. *J. Exp. Med.* **215**(5), 1315-1325 (2018).

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, 09/17/2024

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