

PRODUCT INFORMATION

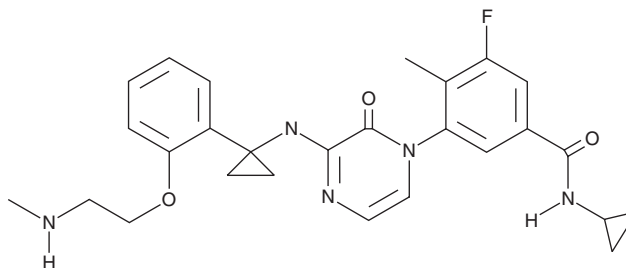


AZD 7624

Item No. 27690

CAS Registry No.: 1095004-78-6
Formal Name: N-cyclopropyl-3-fluoro-4-methyl-5-[3-[[1-[2-[2-(methylamino)ethoxy]phenyl]cyclopropyl]amino]-2-oxo-1(2H)-pyrazinyl]-benzamide

MF: C₂₇H₃₀FN₅O₃
FW: 491.6
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

AZD 7624 is supplied as a solid. A stock solution may be made by dissolving the AZD 7624 in the solvent of choice, which should be purged with an inert gas. AZD 7624 is soluble in acetonitrile and methanol.

Description

AZD 7624 is an inhibitor of p38 α MAPK signaling.¹ AZD 7624 (1 μ M), in combination with the synthetic glucocorticoid dexamethasone (Item No. 11015), prevents LPS-induced increases in the levels of TNF- α and IL-6 secreted from primary human alveolar macrophages isolated from patients with chronic obstructive pulmonary disease (COPD).

Reference

- Cooper, A.E., Luker, T.J., and Schmidt, J.A. Pharmaceutical product comprising a p38 kinase inhibitor and a second active ingredient. *Astrazeneca AB. WO2010/071583A1* (2010).

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, 11/01/2023

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