

PRODUCT INFORMATION

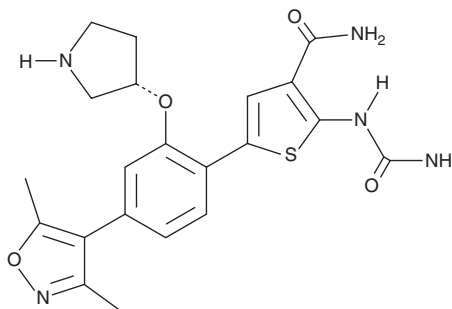


AZD 3264

Item No. 19309

CAS Registry No.: 1609281-86-8
Formal Name: 2-[(aminocarbonyl)amino]-5-[4-(3,5-dimethyl-4-isoxazolyl)-2-[(3S)-3-pyrrolidinyloxy]phenyl]-3-thiophenecarboxamide

MF: C₂₁H₂₃N₅O₄S
FW: 441.5
Purity: ≥98%
UV/Vis.: λ_{max}: 243, 335 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

AZD 3264 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 3264 in the solvent of choice, which should be purged with an inert gas. AZD 3264 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AZD 3264 in these solvents is approximately 30 mg/ml.

Description

AZD 3264 is an inhibitor of NF-κB kinase 2 (IKK2).¹

Reference

1. Murugan, A., Bachu, S., Manjunatha, S.G., *et al.* Exploiting the differential reactivities of halogen atoms: Development of a scalable route to IKK2 inhibitor AZD3264. *Org. Process Res. Dev.* **18(5)**, 646-651 (2014).

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, 11/18/2022

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