

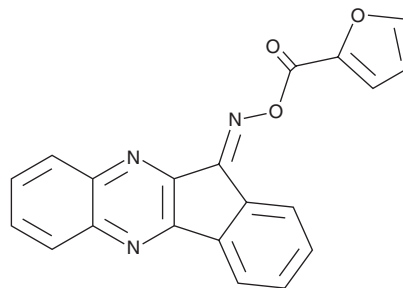
PRODUCT INFORMATION



IQ-3

Item No. 44645

CAS Registry No.: 312538-03-7
Formal Name: O-(2-furanylcarbonyl)oxime
11H-indeno[1,2-b]quinoxalin-11-one
MF: C₂₀H₁₁N₃O₃
FW: 341.3
Purity: ≥98%
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

IQ-3 is supplied as a solid. A stock solution may be made by dissolving the IQ-3 in the solvent of choice, which should be purged with an inert gas. IQ-3 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

IQ-3 is a JNK inhibitor.¹ It is selective for JNKs over a panel of 131 kinases at 10 μM. IQ-3 binds to JNKs (K_d = 0.24, 0.29, and 0.066 μM for JNK1-3), as well as casein kinase 1δ (CK1δ), PI3Kγ, and MAPK-interacting serine/threonine kinase 2 (MKNK2; K_ds = 0.56, 0.43, and 1.2 μM, respectively). It also inhibits NF-κB/AP-1 (IC₅₀ = 1.4 μM in a reporter assay). IQ-3 reduces TNF-α and IL-6 production in human Mono-Mac-6 cells (IC₅₀s = 2.2 and 1.5 μM, respectively) and isolated human peripheral blood mononuclear cells (PBMCs; IC₅₀s = 4.7 and 9.1 μM, respectively), as well as nitric oxide (NO) production in murine J774A.1 macrophages (IC₅₀ = 6.4 μM).

Reference

1. Schepetkin, I.A., Kirpotina, L.N., Khlebnikov, A.I., *et al.* Identification and characterization of a novel class of c-Jun N-terminal kinase inhibitors. *Mol. Pharmacol.* **81(6)**, 832-845 (2012).

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, 12/08/2025

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