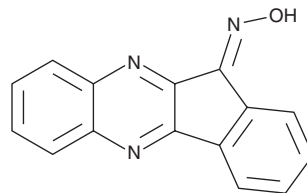


PRODUCT INFORMATION

IQ-1S (free acid)

Item No. 29851

CAS Registry No.: 23146-22-7
Formal Name: 11H-indeno[1,2-b]quinoxalin-11-one, oxime
Synonym: IQ-1
MF: C₁₅H₉N₃O
FW: 247.3
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-1S (free acid) is supplied as a solid. A stock solution may be made by dissolving the IQ-1S (free acid) in the solvent of choice, which should be purged with an inert gas. IQ-1S (free acid) is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of IQ-1S (free acid) in these solvents is approximately 10 and 15 mg/ml, respectively.

IQ-1S (free acid) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IQ-1S (free acid) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. IQ-1S (free acid) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

IQ-1S (free acid) is an inhibitor of NF-κB/AP-1 (IC₅₀ = 2.3 μM in a reporter assay).¹ It inhibits TNF-α and IL-6 production in human Mono-Mac-6 cells (IC₅₀s = 1.3 and 3.8 μM, respectively) and isolated human peripheral blood mononuclear cells (PBMCs; IC₅₀s = 2.6 and 5.6 μM, respectively), as well as nitric oxide (NO) production in murine J774-A.1 macrophages (IC₅₀ = 3.1 μM). IQ-1S (free acid) binds to JNKs (K_d = 0.24, 0.36, and 0.1 μM for JNK1-3), as well as casein kinase 1δ (CK1δ), PI3Kγ, and MAPK-interacting serine/threonine kinase 2 (MKNK2; K_ds = 0.38, 0.47, and 0.92 μM, respectively).

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, 11/03/2022

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