

PRODUCT INFORMATION



BIX02188

Item No. 16238

CAS Registry No.: 334949-59-6
Formal Name: (3Z)-3-[[[3-[(dimethylamino)methyl]phenyl]amino]phenylmethylene]-2,3-dihydro-2-oxo-1H-indole-6-carboxamide

MF: C₂₅H₂₄N₄O₂

FW: 412.5

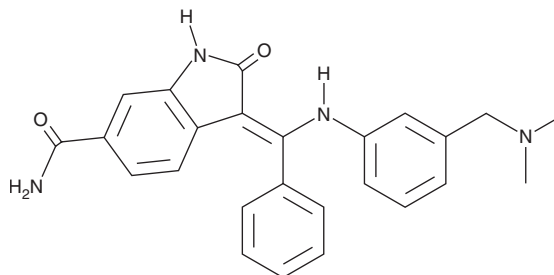
Purity: ≥95%

UV/Vis.: λ_{max}: 210, 285, 384 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

BIX02188 is supplied as a crystalline solid. A stock solution may be made by dissolving the BIX02188 in the solvent of choice, which should be purged with an inert gas. BIX02188 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BIX02188 in these solvents is approximately 20 mg/ml.

BIX02188 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BIX02188 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BIX02188 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BIX02188 is a potent inhibitor of MEK5 (IC₅₀ = 4.3 nM) with reduced activity towards inhibition of ERK5 and TGFβR1 (IC₅₀s = 810 and 1,800 nM, respectively).¹ It does not inhibit the closely related kinases MEK1, MEK2, ERK2, and JNK2 (IC₅₀s = > 6,300 nM).¹ BIX02188 inhibits phosphorylation of ERK5 in a dose-dependent manner without affecting phosphorylation of ERK1/2, p38 MAPK, or JNK1/2.¹

Reference

1. Tataka, R.J., O'Neill, M.M., Kennedy, C.A., *et al.* Identification of pharmacological inhibitors of the MEK5/ERK5 pathway. *Biochem. Biophys. Res. Commun.* **377**(1), 120-125 (2008).

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, 11/30/2022

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