

PRODUCT INFORMATION

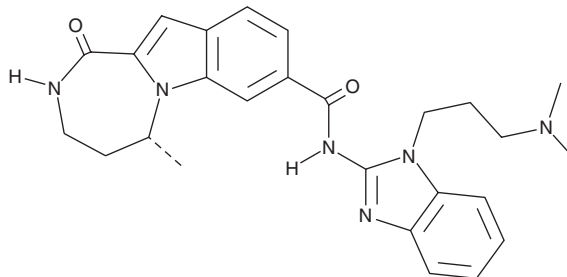


BIX 02565

Item No. 19183

CAS Registry No.: 1311367-27-7
Formal Name: (5R)-N-[1-[3-(dimethylamino)propyl]-1H-benzimidazol-2-yl]-2,3,4,5-tetrahydro-5-methyl-1-oxo-1H-[1,4]diazepino[1,2-a]indole-8-carboxamide

MF: C₂₆H₃₀N₆O₂
FW: 458.6
Purity: ≥98%
UV/Vis.: λ_{max}: 248, 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

BIX 02565 is supplied as a crystalline solid. A stock solution may be made by dissolving the BIX 02565 in the solvent of choice, which should be purged with an inert gas. BIX 02565 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BIX 02565 in these solvents is approximately 1 mg/ml.

BIX 02565 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BIX 02565 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BIX 02565 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

The 90 kDa ribosomal S6 kinases (RSKs) are a group of serine/threonine kinases involved in diverse cellular processes, including growth, survival, and motility. Phosphorylation of the sodium-hydrogen exchanger (NHE) in cardiac tissue by RSK has been shown to activate NHE during cellular stresses, such as myocardial infarction, as a mechanism to maintain pH. BIX 02565 is a RSK2 inhibitor with an IC₅₀ value of 1.1 nM.^{1,2} It also demonstrates off-target binding at multiple adrenergic receptor subtypes that are important for vascular tone and cardiac function (IC₅₀s = 0.052-1.820 μM for adrenergic α_{1A}, α_{1B}, α_{1D}, α_{2A}, β₂, and imidazoline I₂ receptors.^{1,2} BIX 02565 can also inhibit LRRK2 and PRKD1 with IC₅₀ values of 16 and 35 nM, respectively).¹ In rats, BIX 02565 at 1, 3, and 10 mg/kg was shown to significantly decrease heart rate and mean arterial pressure.²

References

1. Kirrane, T.M., Boyer, S.J., Burke, J., *et al.* Indole RSK inhibitors. Part 2: Optimization of cell potency and kinase selectivity. *Bioorg. Med. Chem. Lett.* **22(1)**, 738-742 (2012).
2. Fryer, R.M., Muthukumarana, A., Chen, R.R., *et al.* Mitigation of off-target adrenergic binding and effects on cardiovascular function in the discovery of novel ribosomal S6 kinase 2 inhibitors. *J. Pharmacol. Exp. Ther.* **340(3)**, 492-500 (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, 11/09/2022

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