

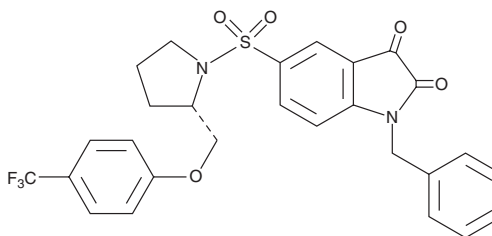
PRODUCT INFORMATION



CAY10406

Item No. 72510

CAS Registry No.: 2108603-03-6
Formal Name: (S)-1-benzyl-5-(1-[2-(phenoxy-*p*-trifluoromethyl)-pyrrolidynyl]-sulfonyl)-isatin
MF: C₂₇H₂₃F₃N₂O₅S
FW: 544.5
Purity: ≥98%
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

CAY10406 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10406 in the solvent of choice. CAY10406 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of CAY10406 in these solvents is approximately .05, 10, and 25 mg/ml, respectively.

CAY10406 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10406 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10406 has a solubility of approximately 0.5 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

CAY10406 is a trifluoromethyl analog of an isatin sulfonamide compound that selectively inhibits caspases 3 and 7.¹ The non-trifluoromethyl compound exhibits K_i values of 1.2 nM and 6 nM for caspases 3 and 7, respectively. For all of the other caspases tested, it is 100 to 1,000 times less potent. Caspases 3 and 7 are 'effector caspases' that are downstream from the initiating steps of apoptosis, and are implicated in the main proteolytic processing of the apoptotic signal. No data is currently available for caspase inhibition by CAY10406.

Reference

1. Lee, D., Long, S.A., Murray, J.H., *et al.* Potent and selective nonpeptide inhibitors of caspases 3 and 7. *J. Med. Chem.* **44**, 2015-2026 (2001).

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, 04/08/2024

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