

PRODUCT INFORMATION

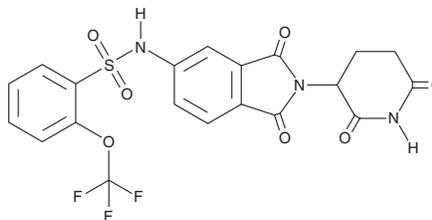


SJ6986

Item No. 34523

CAS Registry No.: 2765625-93-0
Formal Name: N-[2-(2,6-dioxo-3-piperidiny)-2,3-dihydro-1,3-dioxo-1H-isoindol-5-yl]-2-(trifluoromethoxy)benzenesulfonamide

MF: C₂₀H₁₄F₃N₃O₇S
FW: 497.4
Purity: ≥98%
UV/Vis.: λ_{max}: 237 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SJ6986 is supplied as a crystalline solid. A stock solution may be made by dissolving the SJ6986 in the solvent of choice, which should be purged with an inert gas. SJ6986 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SJ6986 in these solvents is approximately 30 mg/ml.

SJ6986 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SJ6986 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SJ6986 has a solubility of approximately 0.20 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SJ6986 is a degrader of G₁ to S phase transition 1 (GSPT1) and GSPT2.¹ It reduces the levels of GSPT1 and GSPT2 in MV4-11 cells in a concentration-dependent manner. SJ6986 inhibits proliferation of MV4-11, MHH-CALL-4, MB002, MB004, and HD-MB03 cells (EC₅₀s = 1.5, 0.4, 726, 336, and 3,583 nM, respectively) and reduces the viability of MV4-11 cells endogenously expressing cereblon but not MV4-11 cereblon knockout cells.

Reference

1. Nishiguchi, G., Keramatnia, F., Min, J., *et al.* Identification of potent, selective, and orally bioavailable small-molecule GSPT1/2 degraders from a focused library of cereblon modulators. *J. Med. Chem.* **64**(11), 7296-7311 (2021).

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, 9/01/2022

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