

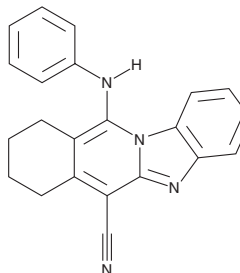
PRODUCT INFORMATION



Usp22i-S02 Item No. 42130

CAS Registry No.: 309735-96-4
Formal Name: 7,8,9,10-tetrahydro-11-(phenylamino)-benzimidazo[1,2-b]isoquinoline-6-carbonitrile
Synonyms: Usp22-IN-1, Usp22 Inhibitor 1, Ubiquitin-specific Protease 22-IN-1, Ubiquitin-specific Protease 22 Inhibitor 1, Ubiquitin-specific Protease 22i-S02

MF: C₂₂H₁₈N₄
FW: 338.4
Purity: ≥95%
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

Usp22i-S02 is supplied as a solid. A stock solution may be made by dissolving the Usp22i-S02 in the solvent of choice, which should be purged with an inert gas. Usp22i-S02 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Usp22i-S02 is an inhibitor of ubiquitin-specific protease 22 (USP22).¹ It increases the levels of ubiquitinated forkhead box protein P3 (Foxp3), a target of Usp22, in primary mouse regulatory T cells (Tregs) when used at concentrations of 10, 15, or 20 μM. *In vivo*, Usp22i-S02 (20 mg/kg per day for five days) reduces tumor volume and increases intratumoral levels of CD4⁺CD8⁺ T cells in an LLC1 murine lung cancer model.

Reference

1. Montauti, E., Weinberg, S.E., Chu, P., *et al.* A deubiquitination module essential for T_{reg} fitness in the tumor microenvironment. *Sci. Adv.* **8(47)**, eabo4116 (2022).

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, 11/11/2024

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