

PRODUCT INFORMATION

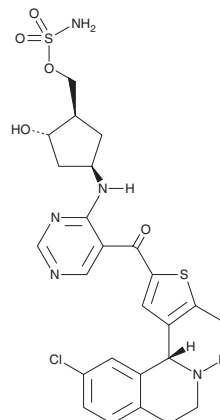


TAK-981

Item No. 32741

CAS Registry No.: 1858276-04-6
Formal Name: [(1R,2S,4R)-4-[[5-[[4-[(1R)-7-chloro-1,2,3,4-tetrahydro-1-isoquinolinyl]-5-methyl-2-thienyl]carbonyl]-4-pyrimidinyl]amino]-2-hydroxycyclopentyl]sulfamic acid, methyl ester

MF: C₂₅H₂₈ClN₅O₅S₂
FW: 578.1
Purity: ≥98%
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

TAK-981 is supplied as a solid. A stock solution may be made by dissolving the TAK-981 in the solvent of choice, which should be purged with an inert gas. TAK-981 is soluble in DMSO.

Description

TAK-981 is an inhibitor of SUMO-activating enzyme (SAE; IC₅₀ = <10 nM).¹ It increases macrophage phagocytosis and natural killer (NK) cell-mediated cytotoxicity of cancer cells *in vitro*.² TAK-981 upregulates the expression of IFN-regulated genes in the blood and tumor tissue in an A20 syngeneic lymphoma mouse model.³ It also decreases intratumor levels of SUMOylated proteins and reduces tumor growth in the same model.

References

1. Duffey, M.O., England, D., Freeze, S., *et al.* Heteroaryl compounds useful as inhibitors of sumo activating enzyme. **WO2016/004136 A1**, (2016).
2. Nakamura, A., Grossman, S., Song, K., *et al.* Abstract 1523: Inhibition of SUMOylation by TAK-981 induces antitumor innate immune responses by modulating macrophage and NK cell function through Type I IFN pathway activation. *Cancer Res.* **79(13)**, 1523 (2019).
3. Berger, A.J., Friedlander, S., Ghasemi, O., *et al.* Abstract 3079: Pharmacodynamic evaluation of the novel SUMOylation inhibitor TAK-981 in a mouse tumor model. *Cancer Res.* **79(13)**, 3079 (2019).

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, 12/13/2022

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