

PRODUCT INFORMATION

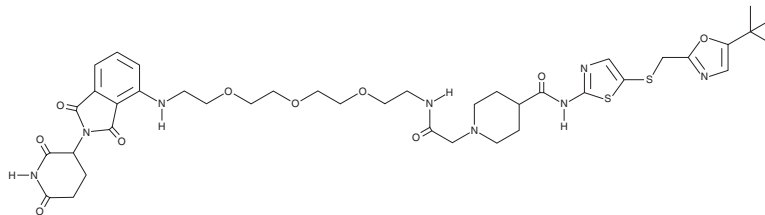


THAL-SNS-032

Item No. 40958

CAS Registry No.: 2139287-33-3
Formal Name: 4-[[[5-[[[5-(1,1-dimethylethyl)-2-oxazolyl]methyl]thio]-2-thiazolyl]amino]carbonyl]-N-[2-[2-[2-[2-[[2-(2,6-dioxo-3-piperidiny)]-2,3-dihydro-1,3-dioxo-1H-isindol-4-yl]amino]ethoxy]ethoxy]ethoxy]ethyl]-1-piperidineacetamide

MF: C₄₀H₅₂N₈O₁₀S₂
FW: 869.0
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

THAL-SNS-032 is supplied as a solid. A stock solution may be made by dissolving the THAL-SNS-032 in the solvent of choice, which should be purged with an inert gas. THAL-SNS-032 is slightly soluble (0.1-1 mg/ml) in methanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

THAL-SNS-032 is a proteolysis-targeting chimera (PROTAC) containing the cyclin-dependent kinase 2 (Cdk2), Cdk7, and Cdk9 inhibitor SNS-032 (Item No. 17904) linked to a cereblon (CRBN) binding moiety.¹ It selectively induces degradation of Cdk9 over Cdk1, Cdk2, and Cdk7 in MOLT-4 cells when used at a concentration of 250 nM. THAL-SNS-032 inhibits human cytomegalovirus (HCMV) replication in infected human foreskin fibroblasts (HFFs; EC₅₀ = 0.025 μM).² It also inhibits viral replication in murine CMV-infected mouse embryonic fibroblasts (MEFs; EC₅₀ = 0.22 μM).

References

1. Olson, C.M., Jiang, B., Erb, M.A., *et al.* Pharmacological perturbation of CDK9 using selective CDK9 inhibition or degradation. *Nat. Chem. Biol.* **14**(2), 163-170 (2018).
2. Hahn, F., Hamilton, S.T., Wangen, C., *et al.* Development of a PROTAC-based targeting strategy provides a mechanistically unique mode of anti-cytomegalovirus activity. *Int. J. Mol. Sci.* **22**(23), 12858 (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, 06/19/2024

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