

PRODUCT INFORMATION



GMP IWR-1-endo Ancillary Material

Item No. 44886

CAS Registry No.: 1127442-82-3

Formal Name: 4-[(3aR,4S,7R,7aS)-1,3,3a,4,7,7a-hexahydro-1,3-dioxo-4,7-methano-2H-isoindol-2-yl]-N-8-quinolinylbenzamide

MF: $C_{25}H_{19}N_3O_3$

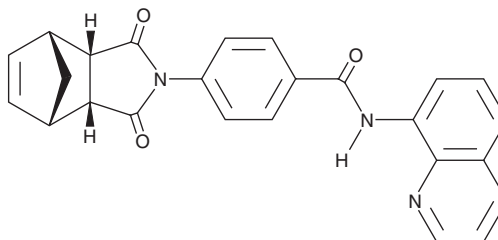
FW: 409.4

Purity: $\geq 98\%$

Supplied as: An amorphous solid

Storage: Room temperature

Stability: **batch specific** Please see the Certificate of Analysis for retest dating.



Description

IWR-1-endo is an inhibitor of the Wnt signaling pathway that acts *via* inhibition of tankyrases and stabilization of axin.^{1,2} It has been used to maintain proliferation of hematopoietic stem cells, promote pluripotency of human embryonic stem cells, and induce differentiation of progenitor cells.³⁻⁵ GMP IWR-1-endo Ancillary Material is intended for use in cell, gene, and tissue engineering applications.

This product has been manufactured and tested in accordance with GMP principles set forth by the ICH and FDA and meets the expectations for commercial cell, gene, and tissue-engineered products in accordance with USP <1043> "Ancillary Materials for Cell, Gene, and Tissue-engineered Products".⁶

References

1. Chen, B., Dodge, M. E., Tang, W., *et al.* Small molecule-mediated disruption of Wnt-dependent signaling in tissue regeneration and cancer. *Nat. Chem. Biol.* **5**(2), 100-107 (2009).
2. Huang, S. M. A., Mishina, Y. M., Liu, S., *et al.* Tankyrase inhibition stabilizes axin and antagonizes Wnt signalling. *Nature* **461**, 614-619 (2009).
3. Kimura, K., Yamamori, S., Hazawa, M., *et al.* Inhibition of canonical Wnt signaling promotes ex vivo maintenance and proliferation of hematopoietic stem cells in zebrafish. *Stem Cells* **40**(9), 831-842 (2022).
4. Choi, H., Oh, D., Kim, M., *et al.* Establishment of porcine embryonic stem cells in simplified serum free media and feeder free expansion. *Stem Cell Res. Ther.* **15**(1), 245 (2024).
5. Ito, A., Ye, K., Onda, M., *et al.* Efficient and robust induction of retinal pigment epithelium cells by tankyrase inhibition regardless of the differentiation propensity of human induced pluripotent stem cells. *Biochem. Biophys. Res. Commun.* **552**, 66-72 (2021).
6. **United States Pharmacopeia.** <1043> Ancillary Materials for Cell, Gene, and Tissue-Engineered Products. In: *USP-NF*. Rockville, MD: United States Pharmacopeial Convention.

WARNING

Ancillary material intended for further processing in accordance with USP <1043>

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, 06/23/2026

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