

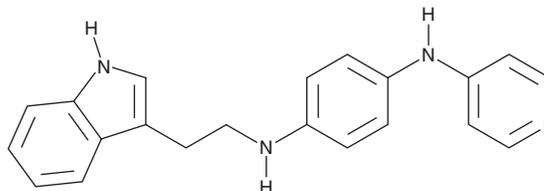
# PRODUCT INFORMATION



**JNJ-26854165**

Item No. 16259

**CAS Registry No.:** 881202-45-5  
**Formal Name:** N<sup>1</sup>-[2-(1H-indol-3-yl)ethyl]-N<sup>4</sup>-4-pyridinyl-1,4-benzenediamine  
**Synonym:** Serdemetan  
**MF:** C<sub>21</sub>H<sub>20</sub>N<sub>4</sub>  
**FW:** 328.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 284 nm  
**Supplied as:** A crystalline 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

JNJ-26854165 is supplied as a crystalline solid. A stock solution may be made by dissolving the JNJ-26854165 in the solvent of choice, which should be purged with an inert gas. JNJ-26854165 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of JNJ-26854165 in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

JNJ-26854165 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JNJ-26854165 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JNJ-26854165 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

The E3 ubiquitin ligase known as murine double minute 2 (MDM2; HDM2 in humans) binds to and then ubiquitinates several proteins, most notably the tumor suppressor p53.<sup>1,2</sup> JNJ-26854165 is an antagonist of MDM2 that suppresses the growth of cancer cell lines expressing wild-type p53 (IC<sub>50</sub> values range from 240-440 nM).<sup>3</sup> It induces p53-mediated transcription culminating in apoptotic death of acute leukemia cells.<sup>3</sup> JNJ-26854165 is orally bioavailable and has anti-proliferative as well as apoptotic actions in various tumor models.<sup>2</sup> As MDM2 interacts with and modulates several targets in addition to p53, JNJ-26854165 affects multiple signaling pathways.<sup>2,4,5</sup>

## References

1. Fulda, S., Rajalingam, K., and Dikic, I. Ubiquitylation in immune disorders and cancer: From molecular mechanisms to therapeutic implications. *EMBO Mol. Med.* **4(7)**, 545-556 (2012).
2. Yuan, Y., Liao, Y.-M., Hsueh, C.-T., et al. Novel targeted therapeutics: Inhibitors of MDM2, ALK and PARP. *J. Hematol. Oncol.* **4(16)**, 1-14 (2011).
3. Kojima, K., Burks, J.K., Arts, J., et al. The novel tryptamine derivative JNJ-26854165 induces wild-type p53- and E2F1-mediated apoptosis in acute myeloid and lymphoid leukemias. *Mol. Cancer Ther.* **9(9)**, 2545-2557 (2010).
4. Lehman, J.A., Hauck, P.M., Gendron, J.M., et al. Serdemetan antagonizes the Mdm2-HIF1a axis leading to decreased levels of glycolytic enzymes. *PLoS One* **8(9)**, 1-5 (2013).
5. Jones, R.J., Gu, D., Bjorklund, C.C., et al. The novel anticancer agent JNJ-26854165 induces cell death through inhibition of cholesterol transport and degradation of ABCA1. *J. Pharmacol. Exp. Ther.* **346(3)**, 381-392 (2013).

### 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, 12/20/2022

## 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