

# Product Information



## (-)-Nutlin-3

Item No. 18585

Sold under license from Hoffman-La Roche

**Formal Name:** (-)-4-(4,5-bis(4-chlorophenyl)-2-(2-isopropoxy-4-methoxyphenyl)-4,5-dihydro-1H-imidazole-1-carbonyl)piperazin-2-one

**Synonym:** Nutlin 3a

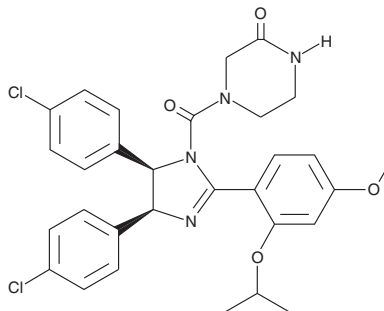
**MF:** C<sub>30</sub>H<sub>30</sub>Cl<sub>2</sub>N<sub>4</sub>O<sub>4</sub>

**FW:** 581.5

**Purity:** ≥98%

**Stability:** ≥2 years at -20°C

**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that (-)-nutlin-3 be stored as supplied at -20°C. It should be stable for at least two years.

(-)-Nutlin-3 is supplied as a crystalline solid. A stock solution may be made by dissolving the (-)-nutlin-3 in an organic solvent purged with an inert gas. (-)-Nutlin-3 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (-)-nutlin-3 in these solvents is approximately 25 mg/ml.

(-)-Nutlin-3 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (-)-nutlin-3 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. (-)-Nutlin-3 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The protein p53, often called the 'guardian of the genome,' is a transcription factor that is activated in response to cellular stress (low oxygen levels, heat shock, DNA damage, etc.) and acts to prevent further proliferation of the stressed cell by promoting cell cycle arrest or apoptosis.<sup>1,2</sup> Its role as a tumor suppressor is evident by the observation that approximately 50% of human tumors have mutated or non-functional p53. Mdm2, a key negative regulator of p53, which is over-expressed in many human tumors, functions by binding to and targeting p53 for proteasomal degradation. Nutlin-3 is a potent inhibitor of p53-Mdm2 interaction.<sup>3</sup> (-)-Nutlin-3 is arbitrarily referred to as enantiomer because it appears as the first peak from chiral purification of racemic nutlin-3 and its absolute stereocenter assignment is not known. It potently inhibits Mdm2-p53 binding with an IC<sub>50</sub> value of 0.09 μM, induces the expression of p53-regulated genes, and exhibits potent antiproliferative activity in cells with functional p53, but not in cells with mutated p53.<sup>3</sup> (-)-Nutlin-3 inhibits the proliferation of exponentially growing human skin and murine fibroblasts with IC<sub>50</sub> values of 2.2 and 1.3 μM, respectively<sup>3</sup>, and induces substantial tumor shrinkage in mice expressing LnCaP, 22Rv1 or MDM cancer cell lines when treated orally with a 200 mg/kg dose twice daily for two weeks.<sup>4</sup>

### References

1. El-Deiry, W.S. The p53 pathway and cancer therapy. *Cancer Journal* **11**, 229-236 (1998).
2. Lane, D.P. and Hupp, T.R. Drug discovery and p53. *Drug Discovery Today* **8(8)**, 347-355 (2003).
3. Vassilev, L.T., Vu, B.T., Graves, B., *et al.* *In vivo* activation of the p53 pathway by small-molecule antagonists of Mdm2. *Science* **303**, 844-848 (2004).
4. Tovar, C., Rosinski, J., Filipovic, Z., *et al.* Small-molecule Mdm2 antagonists reveal aberrant p53 signaling in cancer: Implications for therapy. *Proc. Natl. Acad. Sci. USA* **103(6)**, 1888-1893 (2006).

### Related Products

Nutlin-3 - Item No. 10004372 • (+)-Nutlin-3 - Item No. 10009816

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 01/17/2011

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com