

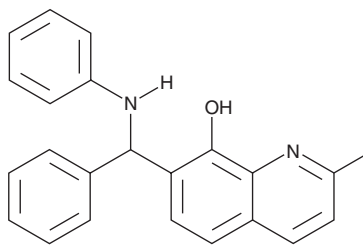
# PRODUCT INFORMATION



## NSC 66811

Item No. 16431

**CAS Registry No.:** 6964-62-1  
**Formal Name:** 2-methyl-7-[phenyl(phenylamino)methyl]-8-quinolinol  
**MF:** C<sub>23</sub>H<sub>20</sub>N<sub>2</sub>O  
**FW:** 340.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 251 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

NSC 66811 is supplied as a crystalline solid. A stock solution may be made by dissolving the NSC 66811 in the solvent of choice, which should be purged with an inert gas. NSC 66811 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NSC 66811 in these solvents is approximately 1, 25, and 33 mg/ml, respectively.

NSC 66811 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NSC 66811 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NSC 66811 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

NSC 66811 is a potent inhibitor of murine double minute 2 (Mdm2) interaction with p53 by binding Mdm2 (K<sub>i</sub> = 120 nM), in this way activating p53.<sup>1</sup> NSC 66811 induces the accumulation of p21, p53, and Mdm2 in human colon cancer cells *in vitro*. Inhibitors of the Mdm2-p53 interaction, including NSC 66811, induce p53- and p21-dependent cell cycle arrest and p53-dependent cell death in tumor cell lines.<sup>2</sup>

### References

1. Lu, Y., Nikolovska-Coleska, Z., Fang, X., *et al.* Discovery of a nanomolar inhibitor of the human murine double minute 2 (MDM2)-p53 interaction through an integrated, virtual database screening strategy. *J. Med. Chem.* **49**(13), 3759-3762 (2006).
2. Shangary, S. and Wang, S. Small-molecule inhibitors of the MDM2-p53 protein-protein interaction to reactivate p53 function: A novel approach for cancer therapy. *Annu. Rev. Pharmacol. Toxicol.* **49**, 223-241 (2016).

#### 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, 10/07/2022

#### 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