

NSC-66811

2-Methyl-7-[Phenyl(phenylamino)methyl]-8-quinolinol

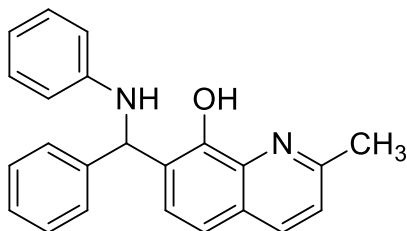
Cat. # SI9850

Background: A cell-permeable, non-peptide inhibitor of MDM2-p53 (HDM2-p53) interactions ($K_i=120\text{nM}$). Treatment of cancer cells activates p53 and accumulation of p21, p53, and MDM2 (HDM2). The structure mimics three p53 residues involved in binding to MDM2 (HDM2).

Application: Cell-based studies of MDM2 (HDM2) and p53 function

Product Information:

CAS No.	6964-62-1
Purity:	>97% (TLC, HPLC), NMR (conforms)
Molecular Weight:	340.4
Physical State:	White Solid
Solubility:	DMSO (25 mg/ml)
Storage:	Store desiccated as supplied at room temperature for 2 years. Store solutions at -20°C for up to 3 months.



Formula: $\text{C}_{23}\text{H}_{20}\text{N}_2\text{O}$

References

1. Lu Y. *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.* 2006, **49**:3759

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved