

Nutlin-3

(±)-4-[4,5-Bis(4-chlorophenyl)-2-(2-isopropoxy-4-methoxy-phenyl)-4,5-dihydro-imidazole-1-carbonyl]-piperazin-2-one

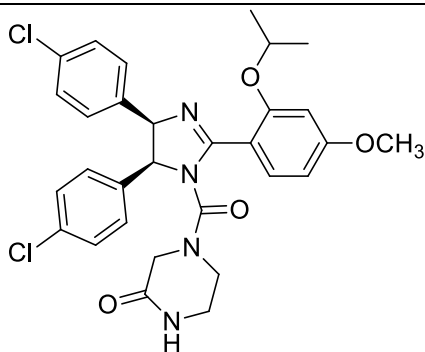
Cat. # SI9870

Background:

Cis-imidazoline analog that inhibits the interaction between p53 and its negative regulator MDM2/HDM2 ($IC_{50} = 0.09 \mu\text{M}$). It induces the expression of p53-regulated genes and exhibits potent antiproliferative activity in cells with functional p53, but not in cells with mutated p53. Nutlin-3 also inhibits the growth of human tumor xenografts in nude mice by 90% at a dose of 200 mg/kg.

Product Information:

CAS No.	548472-68-0
Purity:	>98% (TLC); NMR (Conforms)
Molecular Weight:	581.5
Physical State:	White crystalline solid
Quantity:	5 mg
Solubility:	May be dissolved in DMSO (50 mg/ml)
Storage:	Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 1 month.



Formula: $C_{30}H_{30}Cl_2N_4O_4$

References

1. Vassilev L.T. *et al.* In vivo activation of the p53 pathway by small-molecule antagonists of MDM2. *Science* 2004 **303**(5659):844

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