

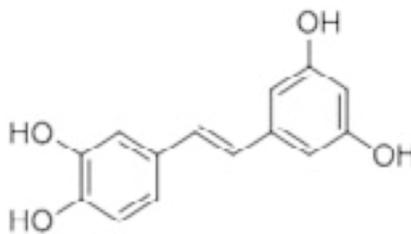
Piceatannol

Catalog Number P009-5MG

Catalog Number P009-25MG

FEATURES

- Activator of SIRT1
- Inhibits NF-κB activation
- Induces apoptosis



ARBOR
ASSAYS

INTRODUCTION

Activator of human deacetylase SIRT1. Anti-inflammatory, immunomodulatory and antiproliferative agent that inhibits p56lck and syk protein tyrosine kinases and inhibits TNF-induced NF-κB activation and gene expression. Piceatannol is a resveratrol analog formed by the cytochrome P450-catalyzed hydroxylation of resveratrol. Piceatannol exhibits potent anticancer properties by inducing apoptosis in BJAB Burkitt-like lymphoma cells with an ED₅₀ value of 25 μM. Piceatannol also exhibits anti-proliferative and anti-inflammatory effects by inhibiting the activity of a wide range of tyrosine and serine/threonine protein kinases and suppressing NF-κB activation through inhibition of IκBα kinase.

FORM: Light pale pink powder

MOLECULAR WEIGHT: 244.25

STORAGE: 4°C

FORMULA: C₁₄H₁₂O₄

CAS NUMBER: 10083-24-6

OTHER NAMES: 4-[(1E)-2-(3,5-Dihydroxyphenyl)ethenyl]-1,2-benzenediol;
3,4,3',5'-Tetrahydroxy-trans-stilbene; Astringenin

USES: Soluble to 10 mg/mL in DMSO or ethanol

REFERENCES:

Wieder, T., Prokop, A., Bagci, B., et al. Piceatannol, a hydroxylated analog of the chemopreventive agent resveratrol, is a potent inducer of apoptosis in the lymphoma cell line BJAB and in primary, leukemic lymphoblasts. *Leukemia* 15, 1735-1742 (2001).

Geahlen, R.L., and McLaughlin, J.L. Piceatannol (3,4,3',5'-tetrahydroxy-trans-stilbene) is a naturally occurring protein-tyrosine kinase inhibitor. *Biochem. Biophys. Res. Comm.* 165:1, 241-245 (1989).

Potter, G.A., Patterson, L.H., Wanogho, E., et al. The cancer preventative agent resveratrol is converted to the anticancer agent piceatannol by the cytochrome P450 enzyme CYP 1B1. *Br. J. Cancer* 86, 774-778 (2002).