Piceatannol



Catalog No: 14036, 14037

Format: 25 mg, 5 mg

Chemical Properties:

MW = 244.2 $C_{14}H_{12}O_4$

CAS 10083-24-6

Physical Properties: Light pink powder **Names:** 3,3',4,5'-Tetrahydroxy-*trans*-stilbene

Pharmacology: A naturally occurring resveratrol analog. Inhibits non-receptor

kinases Syk and Lyk (IC_{50} = ca. 10 M) (ref 1). Stimulates Sirt1 (ref 2).

Solubilization: May be dissolved in DMSO (10 mg/ml); Ethanol (10 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. CJ Seow et al. Eur. J. Pharmacol. 2002, 443:189

2. KT Howitz et al. Nature 2003, 425:191

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 2 months. This product is guaranteed for 6 months from date of arrival.