

Triacetyl resveratrol

Catalog No: 14116, 14117

Format: 250 mg, 50 mg

Chemical Properties:

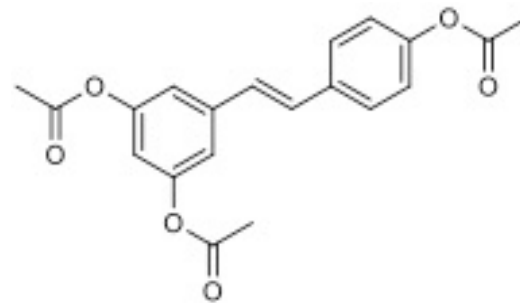
MW = 354.4

C₂₀H₂₈O₆

CAS 42206-94-0

Physical Properties: White powder

Names: Triacetyl resveratrol, 5-[(1E)-2-[4-(acetyloxy)phenyl]ethenyl]-1,3-Benzenediol-1,3-diacetate; 3.5.4'-Tri-O-acetylresveratrol



Pharmacology: A resveratrol prodrug. One method for increasing the half-life of resveratrol *in vivo* is acetylation of the phenolic OH groups. Deacetylation takes place *in vivo* or in intact cells via the action of intracellular esterases releasing active resveratrol.

Solubilization: May be dissolved in DMSO (25 mg/ml); or ethanol (8 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. TC Hsieh *et al. Carcinogenesis* 2011, 32:93
2. TC Hsieh *et al. Int. J. Cancer* 2011, 129:2732

Storage and Guarantee: Store desiccated as supplied at -20° for up to 2 years. Store solutions at -20°C for up to 1 month. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.