Resveratrol



Catalog No: 14022, 14023 Format: 500 mg, 100 mg

Chemical Properties:

MW = 228.2 C₁₄H₁₂O₃ CAS 501-36-0 Physical Properties: Off white powder, MP = 253-255°C **Names:** Resveratrol, 3,4',5-Trihydroxy-trans-stilbene; 5-[(1E)-2-(4-Hydroxyphenyl)ethenyl]-1,3,benzenediol

Pharmacology: Naturally occurring phenolic antioxidant. Displays anti-tumor and anti-platelet effects (ref 1). Activates SIRT1 and increases longevity in yeast (ref 2). Ameliorates ageing-related phenotypes by inhibition of cAMP phosphodiesterases (ref 3).

Solubilization: May be dissolved in DMSO (25 mg/ml); Ethanol (20 mg/ml)

Fluorescent Properties: N/A

Quality Control: >98% (TLC); NMR (Conforms)

References: 1. L Fremont *Life Sci.* 2000, 66:663 2. KT Howitz *et al. Nature* 2003, 425:191

3. S-J Park *et al. Cell* 2012, 148:421

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.

