

PRODUCT DATA SHEET



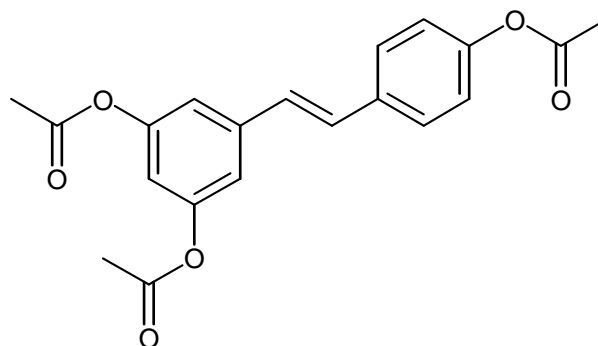
PRODUCT: Triacetyl Resveratrol

CAS NO.: 42206-94-0

CATALOG NO.: FR-119

LOT NO.: Y1069

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 110-112°C)

MOLECULAR FORMULA: C₂₀H₂₈O₆

MOLECULAR WEIGHT: 354.4

PURITY: 98% (TLC: 100% Methylene chloride; R_f = 0.26)

SOLUBILITY: Soluble in; DMSO (25 mg/mL); Ethanol (8 mg/mL warm)

APPLICATION NOTES: A resveratrol prodrug¹. One approach to increasing the half life of resveratrol *in vivo* is acetylation of the phenolic OH groups. Deacetylation presumably takes place *in vivo* or in intact cells via the action of intracellular esterases.

REFERENCES: 1. P. Torres *et al.* *J. Agric. Food Chem.* 2010 **58** 807

STORAGE: Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.

The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.