

PRODUCT DATA SHEET

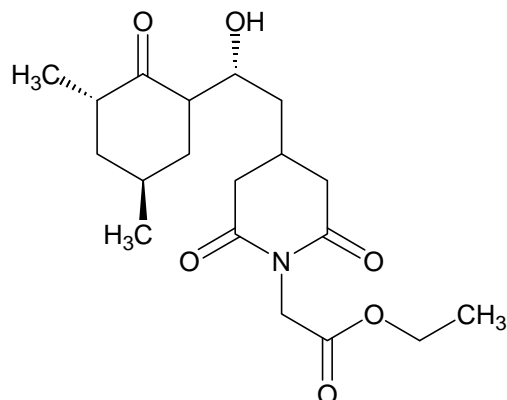


PRODUCT: Cycloheximide-N-ethylethanoate

CATALOG NO.: IM-100

LOT NO.: temp

STRUCTURE:



PHYSICAL APPEARANCE: White solid (m.p. 97-98°C)

MOLECULAR FORMULA: C₁₉H₂₉NO₆

MOLECULAR WEIGHT: 367.4

PURITY: >98%(TLC: 2:1 methylene chloride/ethyl acetate; R_f = 0.43)

SOLUBILITY: Soluble in: DMSO (15 mg/ml); ethanol (3 mg/ml)

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

APPLICATION NOTES: FKBP12 inhibitor (IC₅₀ = 4.4 μM) with neuroregenerative properties. Does not inhibit protein synthesis.

REFERENCES: 1. C.Christner *et al. J.Med.Chem.* 1999 **42** 3615

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.