

PRODUCT DATA SHEET

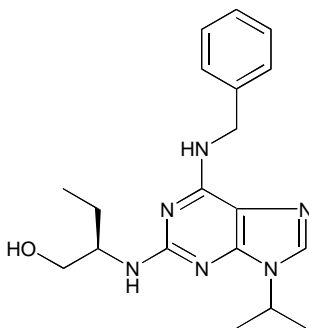


PRODUCT: Roscovitine
2(R)-(1-Ethyl-2-hydroxyethylamino)-
6-benzylamino-9-isopropylpurine

CATALOG NO.: CC-205

LOT#: temp

STRUCTURE:



PHYSICAL APPEARANCE: white solid

MOLECULAR FORMULA: $C_{19}H_{26}N_6O$

MOLECULAR WEIGHT: 354.5

PURITY: 95% (TLC)

SOLUBILITY: Soluble in DMSO

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

APPLICATION NOTES: Cyclin dependent kinase (p34cdc2/cyclin B kinase) inhibitor (IC_{50} = $0.2\ \mu\text{M}$).¹

REFERENCES: 1. L.Havlicek *et al. J.Med.Chem.* 1997 **40** 408

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.