

Roscovitine

Catalog #: 27040

Lot #: 150727

Size: 10 mg

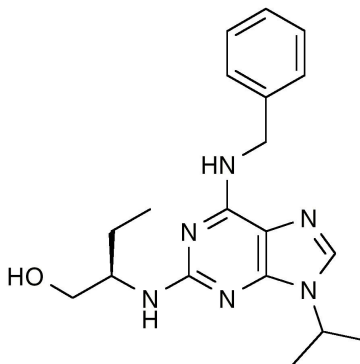
Structure:

CAS Registry #: 186692-46-6

Purity: ≥ 99%

Chemical Formula: C₁₉H₂₆N₆O

Molecular Weight: 354.5 Da



Description: Roscovitine is a potent inhibitor of p34cdc2 (IC₅₀ = 0.2 μM). Also inhibits p33cdk2 and p33cdk5.

Appearance: White to off-white powder.

Solubility: Prepare initial stock solutions in DMSO (soluble at 200 mg/ml) or ethanol (soluble at 200 mg/ml) for in vitro use; solubility in water is low.

Storage/Stability: Store at or below -20°C. Solid form is stable at least 12 months from date of receipt, when stored as directed. Do not store aqueous solutions for more than one day.

Quality Control: The purity was determined by HPLC.

References: Havlicek, L., et al. J. Med. Chem. 40: 408-412 (1997).