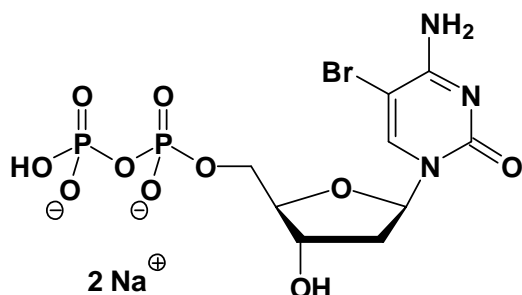


5Br-dCDP

5-Bromo-2'-deoxy-cytidine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-133S	50 Units
NU-133L	250 Units



Cat. No.: NU-133

Molecular Formula: $C_9H_{12}BrN_3O_{10}P_2$ (Anion)

Molecular Weight: 464.05 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Spectroscopic properties: λ_{max} 287 nm; ϵ 6000

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

1 unit = 1 μ l of a 10 mM solution