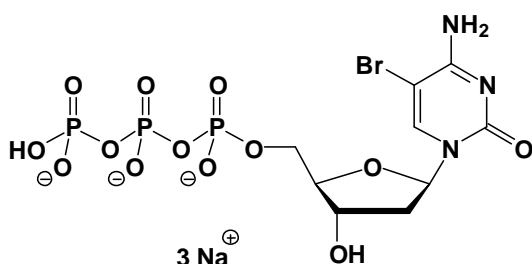


5Br-dCTP

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

Cat. No.	Amount
NU-132S	50 Units
NU-132L	250 Units



Cat. No.: NU-132

Molecular Formula: C₉ H₁₂ BrN₃O₁₃P₃ (Anion)

Molecular Weight: 543.03 (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 μl of a 10 mM solution