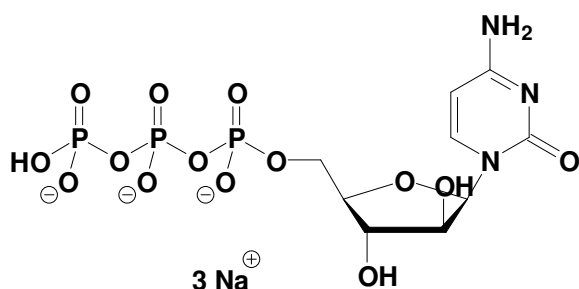


ara-CTP

Cytidine-arabinofuranoside-5'-triphosphate, Sodium salt
(Cytarabine triphosphate)

Cat. No.	Amount
NU-1170S	100 Units
NU-1170L	500 Units



Cat. No.: NU-1170

Molecular Formula: C₉H₁₃N₃O₁₄P₃ (Anion)

Molecular Weight: 480.13 (Anion)

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

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.

* 1 unit = 1 µl of a 10 mM solution

For research use only!