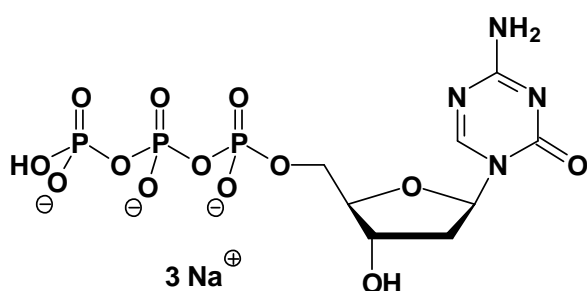


5-Aza-dCTP

5-Aza-2'-deoxy-cytidine-5'-triphosphate, Sodium salt
(Decitabine Triphosphate)

Cat. No.	Amount
NU-1118S	20 Units
NU-1118L	100 Units



Cat. No.: NU-1118

Molecular Formula: C₈H₁₂N₄O₁₃P₃ (Anion)

Molecular Weight: 465.12 (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.

For research use only!

1 unit = 1 µl of a 10 mM solution

Selected References:

Parker *et al.* (1987) Inhibition of DNA primase by nucleoside triphosphates and their arabinofuranosyl analogs. *Mol. Pharmacol.* **31** (2):146.

McIntosh *et al.* (1985) Synthesis and characterization of poly[d(G-z5C)]. B-Z transition and inhibition of DNA methylase. *Biochemistry* **24** (18):4806.

Momparler *et al.* (1984) Kinetic interaction of 5-AZA-2'-deoxycytidine-5'-monophosphate and its 5'-triphosphate with deoxycytidylate deaminase. *Mol. Pharmacol.* **25** (3):436.

Bouchard *et al.* (1983) Incorporation of 5-Aza-2'-deoxycytidine-5'-triphosphate into DNA. Interactions with mammalian DNA polymerase alpha and DNA methylase. *Mol. Pharmacol.* **24** (1):109.

Bouchard, J., Momparler, R. L. (1983) *Molecular Pharmacology* **24-1**:109-114

Richard, L. *et al.* (2005) *Anti-Cancer Drugs.* **16-3**: 301-308