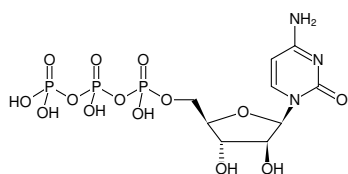




ara-Cytidine-5'-triphosphate (ara-CTP)

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

Cat. No.	Amount
NU-1170S	100 µl (10 mM)
NU-1170L	5 x 100 µl (10 mM)



Structural formula of ara-Cytidine-5'-triphosphate (ara-CTP)

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₉H₁₆N₃O₁₄P₃ (free acid)

Molecular Weight: 483.16 g/mol (free acid)

Exact Mass: 482.98 g/mol (free acid)

CAS#: 13191-15-6

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 10 mM - 11 mM

pH: 7.5 ± 0.5

Spectroscopic Properties: λ_{max} 271 nm, ε 8.9 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)

Applications:

Treatment of acute myeloid leukemia^[1]

Inhibition of DNA synthesis by incorporation^[2]

Detection with isocratic HPLC^[3]

Inhibition of DNA ligases and polymerases^[4]

Selected References:

- [1] Negoro *et al.* (2011) Characterization of cytarabine-resistant leukemic cell lines established from five different blood cell lineages using gene expression and proteomic analyses. *Int. J. Oncol.* **38** (4):911.
- [2] Yamauchi *et al.* (2009) Intracellular cytarabine triphosphate production correlates to deoxycytidine kinase/cytosolic 5'-nucleotidase II expression ratio in primary acute myeloid leukemia cells. *Biochem. Pharmacol.* **77** (12):1780.
- [3] Nishi *et al.* (2006) A new, simple method for quantifying gemcitabine triphosphate in cancer cells using isocratic high-performance liquid chromatography. *Cancer Sci.* **97** (11):1274.
- [4] Colón-Cesario *et al.* (2006) An inhibitor of DNA recombination blocks memory consolidation, but not reconsolidation, in context fear conditioning. *J. Neurosci.* **26** (20):5524.