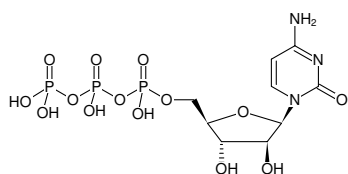




## ara-Cytidine-5'-triphosphate (ara-CTP)

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 research use only!

**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:** C<sub>9</sub>H<sub>16</sub>N<sub>3</sub>O<sub>14</sub>P<sub>3</sub> (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<sub>max</sub> 271 nm, ε 8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.0)

### Applications:

Treatment of acute myeloid leukemia<sup>[1]</sup>

Inhibition of DNA synthesis by incorporation<sup>[2]</sup>

Detection with isocratic HPLC<sup>[3]</sup>

Inhibition of DNA ligases and polymerases<sup>[4]</sup>

### 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.