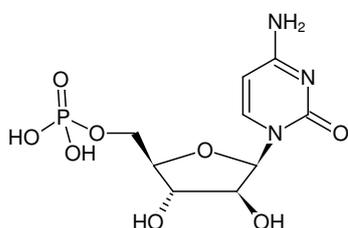



**ara-Cytidine-5'-monophosphate (ara-CMP)**

Cytarabine monophosphate, Sodium salt

Cat. No.	Amount
NU-872	2 mg



Structural formula of ara-Cytidine-5'-monophosphate (ara-CMP)

**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:** C<sub>9</sub>H<sub>14</sub>N<sub>3</sub>O<sub>8</sub>P (free acid)

**Molecular Weight:** 323.20 g/mol (free acid)

**Exact Mass:** 323.05 g/mol (free acid)

**CAS#:** 7075-11-8

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Selected References:**

 Guo *et al.* (2009) Expression of ABC-type nucleotide exporters in blasts of adult acute myeloid leukemia: relation to long-term survival. *Clin Cancer Res* **15** (5):1762.

 Gouy *et al.* (2006) Quantification of 5'-monophosphate cytosine arabinoside (Ara-CMP) in cell extracts using liquid chromatography-electrospray mass spectrometry. *Analytica Chimica Acta* **566** (2):178.

 Hsu *et al.* (2005) Phosphorylation of cytidine, deoxycytidine, and their analog monophosphates by human UMP/CMP kinase is differentially regulated by ATP and magnesium. *Molecular Pharmacology* **67** (3):806.

 Pasti *et al.* (2003) Reaction of human UMP-CMP kinase with natural and analog substrates. *European Journal of Biochemistry* **270** (8):1784.

 Chiramonte *et al.* (2001) Inhibition of CMP-Sialic Acid Transport into Golgi Vesicles by Nucleoside Monophosphates. *Biochemistry* **40** (47):14260.