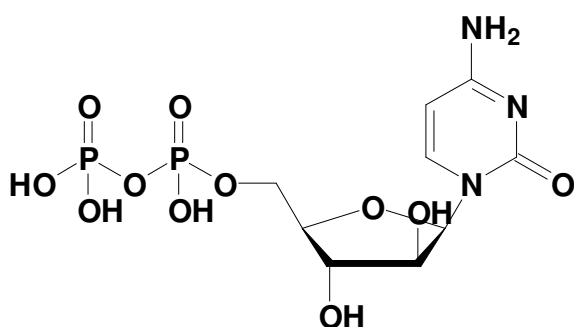


## Ara-cytidine-5'-diphosphate (ara-CDP)

Cytidine-arabinofuranoside-5'-diphosphate, Sodium salt

(Cytarabine diphosphate)

Cat. No.	Amount
NU-1171S	50 Units
NU-1171L	250 Units



**Cat. No.:** NU-1171

**Molecular Formula:**  $C_9H_{15}N_3O_{11}P_2$  (free acid)

**Molecular Weight:** 403.17 (free acid)

**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  $\mu$ l of a 10 mM solution

**For research use only!**