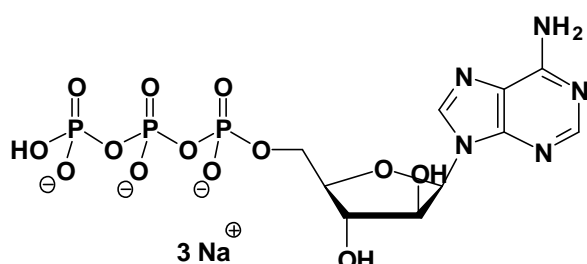


ara-ATP

Adenine-arabinofuranoside-5'-triphosphate, Sodium salt
(Vidarabine triphosphate)

Cat. No.	Amount
NU-1111S	50 Units
NU-1111L	250 Units



Cat. No.: NU-1111

Molecular Formula: C₁₀H₁₃N₅O₁₃P₃ (Anion)

Molecular Weight: 504.16 (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:

Genini *et al.* (2000) Nucleotide requirements for the in vitro activation of the apoptosis protein-activating factor-1-mediated caspase pathway. *J. Biol. Chem.* **275** (1):29.

Ghoshal *et al.* (1991) Ara-ATP impairs 3'-end processing of pre-messenger-RNAs by inhibiting both cleavage and polyadenylation. *Nucleic Acids Res.* **19** (21):5871.

Biron *et al.* (1991) Selective anabolism of 6-methoxypurine arabinoside in Varicella-zoster virus-infected cells. *Antimicrob. Agents Ch.* **35** (10):2116.

Demiranda *et al.* (1991) Anabolic pathway of 6-methoxypurine arabinoside in cells infected with Varicella-zoster virus. *Antimicrob. Agents Ch.* **35** (10):2121.

Plunkett *et al.* (1981) Cellular concentrations of dATP and ara-ATP in peripheral-blood leukemic-cells (PBC) and erythrocytes (RBC) from patients treated with deoxycoformycin (DCF) or Arabinosyladenine (ara-A). *P. Am. Assoc. Canc. Res.* **22** (Mar):177.

Shewach *et al.* (1981) Relationship between ara-ATP concentration, DNA synthetic capacity (DSC) and viability in cho cells. *P. Am. Assoc. Canc. Res.* **22** (Mar):255.

Spater *et al.* (1981) Neuronal phosphatase-activities with ara-AMP and ara-ATP as substrates. *J. Histochem. Cytochem.* **29** (6):693.

Chanda *et al.* (1980) Inhibition of vesicular stomatitis-virus transcriptase invitro by phosphonoformate and ara-ATP. *Virology* **107** (2):562.

Dennis *et al.* (1980) RNA-polymerase - potent linear competitive-inhibition by D Arabinose-5-triphosphate compared to non-inhibition by 5' ara-ATP. *Biochem. Bioph. Res. Co.* **94** (1):205.

Shewach *et al.* (1979) Effect of deoxycoformycin (DCF) on the toxicity and metabolism of 9-beta-d-arabinofuranosyladenine (ara-A) and its 5'-triphosphate (ara-ATP) in cho cells. *P. Am. Assoc. Canc. Res.* **20** (Mar):183.

Chanda *et al.* (1979) Inhibition of invitro RNA-transcription of vesicular stomatitis-virus by phosphonoformate and ara-ATP. *Fed. Proc.* **38** (3):814.

Rose *et al.* (1978) Selective-inhibition of RNA polyadenylation by ara-ATP invitro - possible mechanism for anti-viral action of ara-A. *Biochem. Bioph. Res. Co.* **81** (4):1418.

Ortiz (1972) Inhibition of escherichia-coli adenylyl-cyclase by ara ATP. *Biochem. Bioph. Res. Co.* **46** (4):1728.