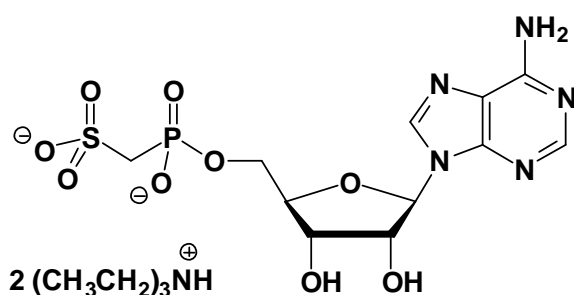


β -methylene-APS

β -Methylene-adenosine-5'-phosphosulfate, Triethylammonium salt

Cat. No.	Amount
NU-701S	50 Units
NU-701L	250 Units



Cat. No.: NU-701

Molecular Formula: C₁₁H₁₄N₅O₉PS (Anion)

Molecular Weight: 423.29 (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 μ l of a 10 mM solution

Selected References:

Menendez-Arias *et al.* (2005) Suppression of Multidrug-resistant HIV-1 Reverse Transcriptase Primer Unblocking Activity by α -Phosphate-modified Thymidine Analogues. *J. Mol. Biol.* **349** (3):451

Kanno *et al.* (1996) Adenosine 5'-phosphosulfate sulfotransferase from the marine macroalga *Porphyra yezoensis* Ueda (Rhodophyta): Stabilization, purification, and properties. *Planta* **198** (3):440.

Lyle *et al.* (1994) Kinetic mechanism of ATP-sulfurylase from rat chondrosarcoma. *Biochem. J.* **301**:349.

Callahan *et al.* (1989) Synthesis and properties of a nonhydrolyzable adenosine phosphosulfate analog. *Analyt. Biochem.* **177** (1):67.