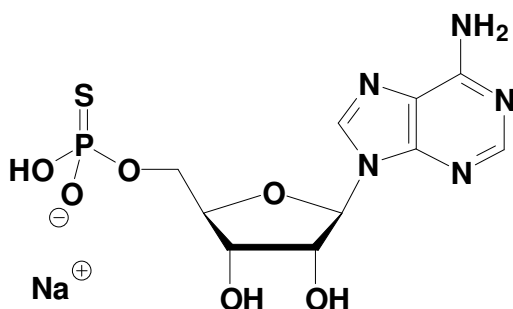


AMP α S

Adenosine-5'-(α -thio)-monophosphate, Sodium salt

Cat. No.	Amount
NU-1161S	100 Units
NU-1161L	500 Units



Cat. No.: NU-1161

Molecular Formula: C₁₀H₁₃N₅O₆PS (Anion)

Molecular Weight: 362.27 (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.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!