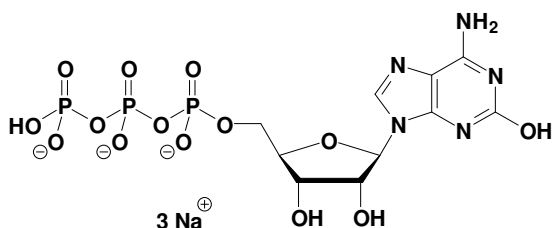


2-Hydroxy-ATP

2-Hydroxy-Adenosine-5'-triphosphate, Sodium salt
(*iso*-GTP)

Cat. No.	Amount
NU-1173S	50 Units
NU-1173L	250 Units



Cat. No.: NU-1173

Molecular Formula: C₁₀H₁₃N₅O₁₄P₃ (Anion)

Molecular Weight: 520.15 (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 µl of a 10 mM solution

For research use only!