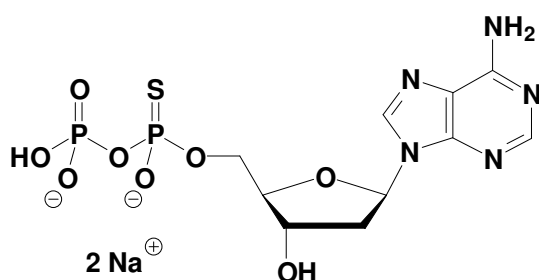


dADP α S

2'-Deoxy-Adenosine-5'-(α -thio)-diphosphate, Sodium salt

Cat. No.	Amount
NU-1165S	20 Units
NU-1165L	100 Units



Cat. No.: NU-1165

Molecular Formula: C₁₀H₁₃N₅O₈P₂S (Anion)

Molecular Weight: 425.24 (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!