

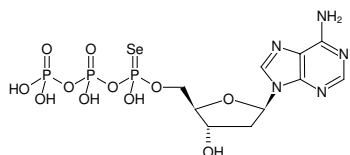
**dATPαSe**

2'-Deoxyadenosine-5'-(α-seleno)-triphosphate, Sodium salt (Mixture of Rp and Sp isomers)

Cat. No.	Amount
NU-267S	10 µl (100 mM)
NU-267L	5 x 10 µl (100 mM)

Selected References:

Hu *et al.* (2019) Synthesis of Selenium-Triphosphates (dNTPαSe) for More Specific DNA Polymerization. *Angew. Chem. Int. Ed.* **58** (23):7835.



Structural formula of dATPαSe

For general laboratory use.**Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₀H₁₆N₅O₁₁P₃Se (free acid)**Molecular Weight:** 554.15 g/mol (free acid)**Exact Mass:** 554.92 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 259 nm, ε 15.4 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)