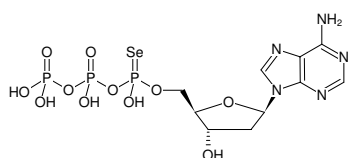


**dATPaSe**2'-Deoxyadenosine-5'-( $\alpha$ -seleno)-triphosphate, Sodium salt (1:1 Mixture of Rp and Sp isomers)

Cat. No.	Amount
NU-267S	10 $\mu$ l (100 mM)
NU-267L	5 x 10 $\mu$ l (100 mM)



Structural formula of dATPaSe

**For in vitro use only!****Shipping:** shipped on blue ice**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:** C<sub>10</sub>H<sub>16</sub>N<sub>5</sub>O<sub>11</sub>P<sub>3</sub>Se (free acid)**Molecular Weight:** 554.15 g/mol (free acid)**Exact Mass:** 554.92 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  259 nm,  $\epsilon$  15.4 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**Nick Translation with Taq Polymerase in-house data**Selected References:**

Hu *et al.* (2019) Synthesis of Selenium<sup>VI</sup> Triphosphates (dNTPaSe) for More Specific DNA Polymerization. *Angew. Chem. Int. Ed.* **58** (23):7835.