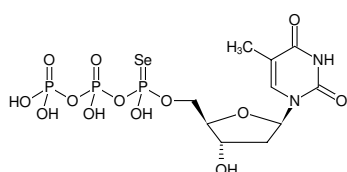


**dTTPαSe**

2'-Deoxythymidine-5'-(α-seleno)-triphosphate, Sodium salt (1:1 Mixture of Rp and Sp isomers)

Cat. No.	Amount
NU-270S	10 µl (100 mM)
NU-270L	5 x 10 µl (100 mM)



Structural formula of dTTPαSe

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:** C₁₀H₁₇N₂O₁₃P₃Se (free acid)**Molecular Weight:** 545.14 g/mol (free acid)**Exact Mass:** 545.91 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 100 mM - 110 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 267 nm, ε 9.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:**Nick Translation with Taq Polymerase in-house data**Selected References:**

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