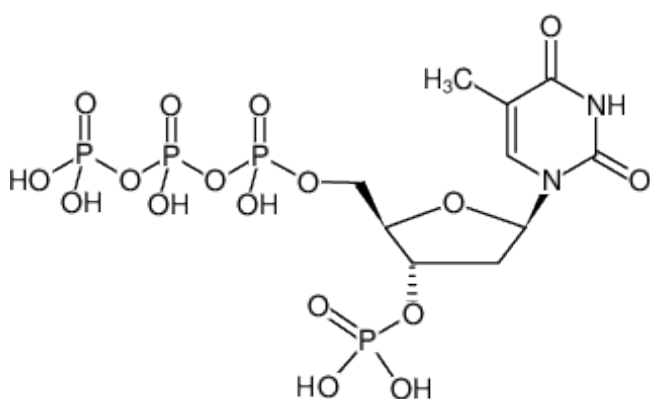


**3'-Phosphate-dTTP**

3'-Phosphate-2'-deoxythymidine-5'-triphosphate, Triethylammonium salt

Cat. No.	Amount
NU-1224S	200 µl (10 mM)
NU-1224L	5 x 200 µl (10 mM)



Structural formula of 3'-Phosphate-dTTP

**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.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>18</sub>N<sub>2</sub>O<sub>17</sub>P<sub>4</sub> (free acid)**Molecular Weight:** 562.15 g/mol (free acid)**Exact Mass:** 561.96 g/mol (free acid)**Purity:** ≥ 98 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ<sub>max</sub> 267 nm, ε 9.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)