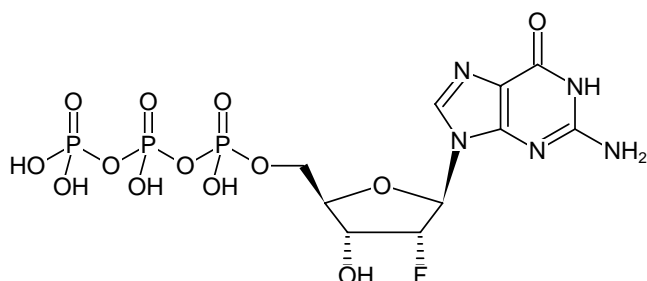


**2'-Fluoro-dGTP**

(2'F-dGTP)

2'-Fluoro-2'-deoxyguanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1216S	50 µl (100 mM)
NU-1216L	5 x 50 µl (100 mM)



Structural formula of 2'-Fluoro-dGTP

For research use only!**Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark and dry

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery**Molecular Formula:** C₁₀H₁₅FN₅O₁₃P₃ (free acid)**Molecular Weight:** 525.17 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} 252 nm; ε 14.2 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.0)