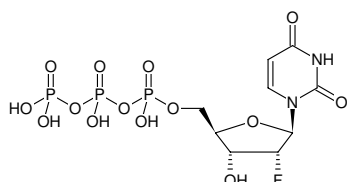


**2'-Fluoro-dUTP**

(2'F-dUTP)

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

Cat. No.	Amount
NU-1215S	50 µl (100 mM)
NU-1215L	5 x 50 µl (100 mM)



Structural formula of 2'-Fluoro-dUTP

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₉H₁₄FN₂O₁₄P₃ (free acid)**Molecular Weight:** 486.13 g/mol (free acid)**Exact Mass:** 485.96 g/mol (free acid)**CAS#:** 66840-02-6**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} 262 nm, ε 10.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)