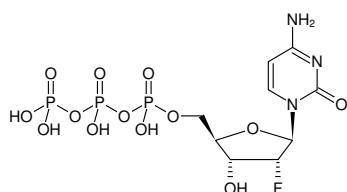


**2'-Fluoro-dCTP**

(2'F-dCTP)

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

Cat. No.	Amount
NU-1214S	50 µl (100 mM)
NU-1214L	5 x 50 µl (100 mM)



Structural formula of 2'-Fluoro-dCTP

**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:** C<sub>9</sub>H<sub>15</sub>FN<sub>3</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 485.15 g/mol (free acid)**Exact Mass:** 484.98 g/mol (free acid)**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<sub>max</sub> 271 nm, ε 9.1 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Related Products:**

HighYield T73M Aptamer Synthesis Kit (2'F-dCTP), #RNT-302

**Selected References:**Ono *et al.* (1997) 2'-Fluoro modified nucleic acids: polymerase-directed synthesis, properties and stability to analysis by matrix-assisted laser desorption/ionization mass spectrometry. *Nucleic Acids Res.* **25 (22)**:4581.