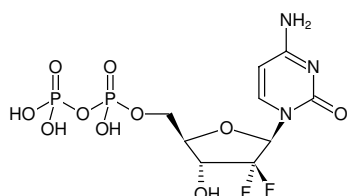


**Gemcitabine-5'-diphosphate**

Sodium Salt

2',2'-Difluorocytidine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-1608S	20 µl (10 mM)
NU-1608L	5 x 20 µl (10 mM)



Structural formula of Gemcitabine-5'-diphosphate

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:** C₉H₁₃N₃O₁₀P₂F₂ (free acid)**Molecular Weight:** 423.16 g/mol (free acid)**Exact Mass:** 423.00 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5**Spectroscopic Properties:** λ_{max} 271 nm, ε 9.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)