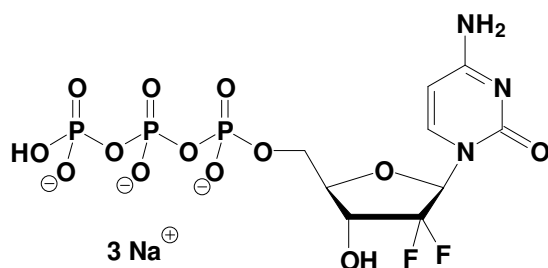


Gemcitabine-Triphosphate

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

Cat. No.	Amount
NU-1607S	50 Units
NU-1607L	250 Units



Cat. No.: NU-1607

Molecular Formula: $C_9H_{11}N_3O_{13}P_3F_2$ (Anion)

Molecular Weight: 500.11 (Anion)

Purity: > 95%, clear aqueous solution, pH 7.5

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

* 1 unit = 1 μ l of a 10 mM solution

For research use only!