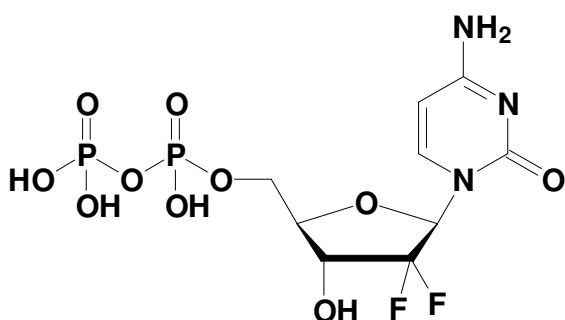


Gemcitabine-Diphosphate

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

Cat. No.	Amount
NU-1608S	20 Units
NU-1608L	100 Units



Cat. No.: NU-1608

Molecular Formula: $C_9H_{13}N_3O_{10}P_2F_2$ (free acid)

Molecular Weight: 423.16 (free acid)

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

Spectroscopic Properties: λ_{max} 271 nm;
 ϵ 9,100 $M^{-1}cm^{-1}$ Tris HCl 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!