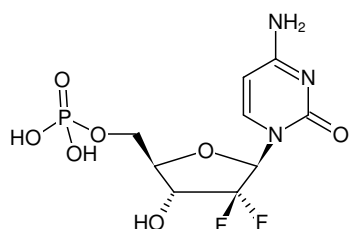




Gemcitabine-5'-monophosphate

Sodium salt

Cat. No.	Amount
NU-871	2 mg



Structural formula of Gemcitabine-5'-monophosphate

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: C₉H₁₂F₂N₃O₇P (free acid)

Molecular Weight: 343.18 g/mol (free acid)

Exact Mass: 343.04 g/mol (free acid)

CAS#: 116371-67-6

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{abs} 271 nm, ε 9.1 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Selected References:

Aksoy *et al.* (2009) Cytosolic 5'-nucleotidase III (NT5C3): gene sequence variation and functional genomics. *Pharmacogenet Genomics* **19** (8):567.

Veltkamp *et al.* (2008) New insights into the pharmacology and cytotoxicity of gemcitabine and 2',2'-difluorodeoxyuridine. *Mol Cancer Ther* **7** (8):2415.

Kocabas *et al.* (2008) Gemcitabine pharmacogenomics: deoxycytidine kinase and cytidylate kinase gene resequencing and functional genomics. *Drug Metab Dispos* **36** (9):1951.