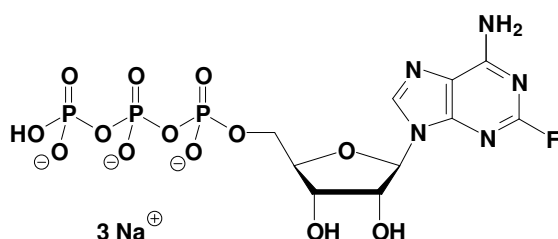


2F-ATP

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

Cat. No.	Amount
NU-145S	20 Units
NU-145L	100 Units



Cat. No.: NU-145

Molecular Formula: $\text{C}_{10}\text{H}_{12}\text{N}_5\text{O}_{13}\text{P}_3\text{F}$ (Anion)

Molecular Weight: 522.15 (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 μl of a 10 mM solution

For research use only!