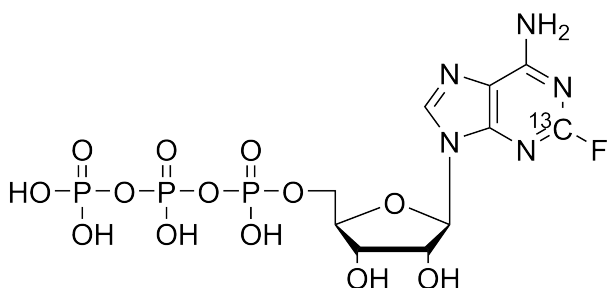


**2-F-2-¹³C-ATP**[2-F, 2-¹³C]-adenosine-5'-triphosphate, lithium salt

Cat. No.	Amount
NU-IT-101-S	100 µl (100 mM)
NU-IT-101-L	5 x 100 µl (100 mM)

Isotopic enrichment: ≥ 98% (Isotopic Enrichment of Precursor)**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:** C₉¹³CH₁₅FN₅O₁₃P₃ (free acid)**Molecular Weight:** 526.16 g/mol (free acid)**Exact Mass:** 525.99 g/mol (free acid)**CAS#:** 1492-62-2 (unlabeled, free acid)**Purity:** ≥ 90 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 100 mM - 110 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 260 nm, ε 13.4 L mmol⁻¹ cm⁻¹ (aqueous solution)