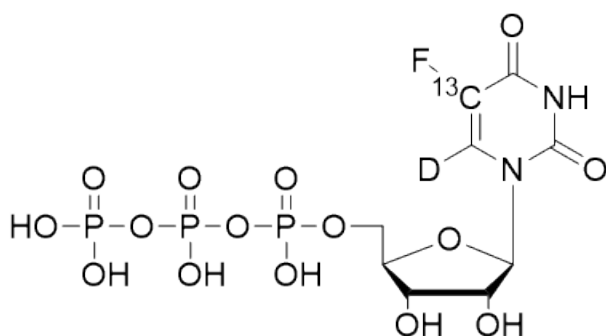


**5-F-5-¹³C-6-D-UTP**[5-F, 5-¹³C, 6-D]-uridine-5'-triphosphate, lithium salt

Cat. No.	Amount
NU-IT-102-S	100 µl (100 mM)
NU-IT-102-L	5x 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:** C₈¹³CH₁₃DFN₂O₁₅P₃ (free acid)**Molecular Weight:** 504.13 g/mol (free acid)**Exact Mass:** 503.97 g/mol (free acid)**CAS#:** 3828-96-4 (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} 270 nm, ε 8.8 L mmol⁻¹ cm⁻¹ (aqueous solution)