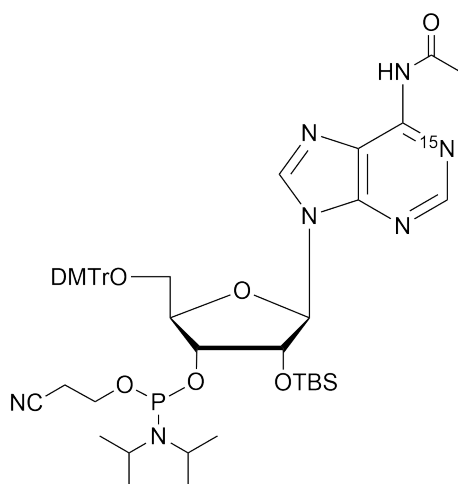


**1-¹⁵N-rA-CEP**[1-¹⁵N]-N6-Acetyl-5'-O-DMT-2'-O-TBDMS-adenosine-3'-O-CE phosphoramidite

Cat. No.	Amount
N-IT-005-S	10 mg
N-IT-005-L	50 mg

Applications:

RNA oligonucleotide synthesis (bottle configured for ABI)

Isotopic enrichment: ≥ 98% (Isotopic Enrichment of Precursor)**For general laboratory use.****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C₄₈H₆₄N₆¹⁵NO₈PSi**Molecular Weight:** 927.13 g/mol**Exact Mass:** 926.43 g/mol**CAS#:** 1220889-84-8 (unlabeled)**Purity:** ≥ 95 % (31P-NMR)**Form:** solid**Color:** white of off-white