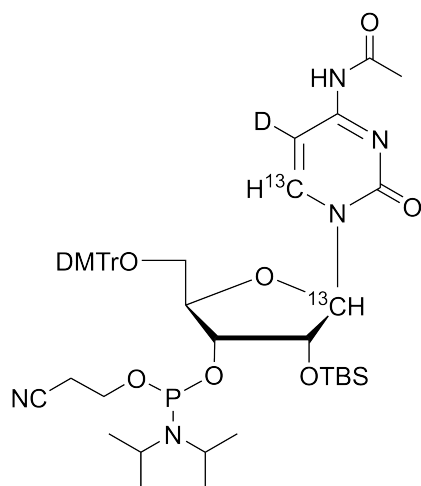


**5-D-1',6-¹³C₂-rC-CEP**[5-D, 1,6-¹³C]-N₄-Acetyl-5'-O-DMT-2'-O-TBDMS-cytidine-3'-O-CE phosphoramidite

Cat. No.	Amount
N-IT-004-S	10 mg
N-IT-004-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:** C₄₅¹³C₂H₆₃DN₅O₉PSi**Molecular Weight:** 905.1 g/mol**Exact Mass:** 904.43 g/mol**CAS#:** 121058-88-6 (unlabeled)**Purity:** ≥ 95 % (31P-NMR)**Form:** solid**Color:** white of off-white