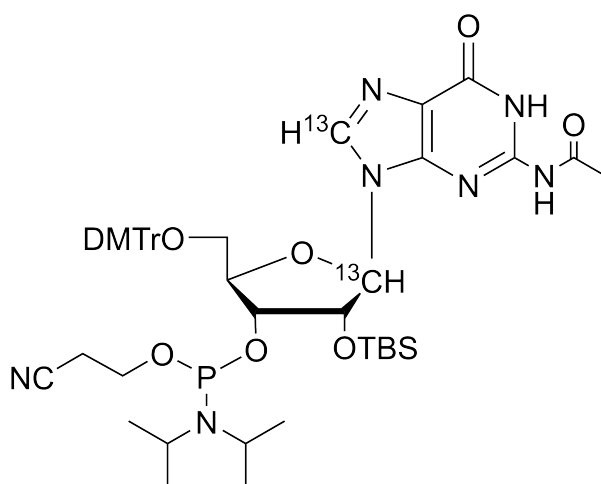


**1',8-<sup>13</sup>C<sub>2</sub>-rG-CEP**[1',8-<sup>13</sup>C]-N<sup>2</sup>-Acetyl-5'-O-DMT-2'-O-TBDMS-guanosine-3'-O-CE phosphoramidite

Cat. No.	Amount
N-IT-002-S	10 mg
N-IT-002-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:** C<sub>46</sub><sup>13</sup>C<sub>2</sub>H<sub>64</sub>N<sub>7</sub>O<sub>9</sub>PSi**Molecular Weight:** 944.12 g/mol**Exact Mass:** 943.43 g/mol**CAS#:** 121058-88-6 (unlabeled)**Purity:** ≥ 95 % (31P-NMR)**Form:** solid**Color:** white of off-white