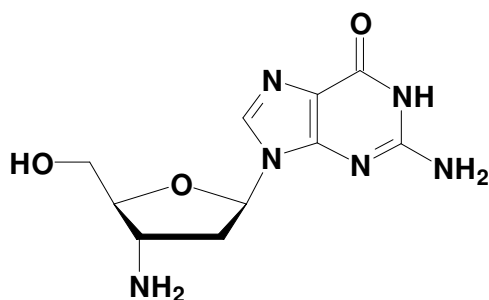


## 3'-Amino-2',3'-Dideoxyguanosine

9-(3-Amino-2,3-dideoxy- $\beta$ -D-ribofuranosyl)-guanine

| Cat. No. | Amount |
|----------|--------|
| N-1010   | 1 g    |



**Cat. No.:** N-1000

**Molecular Formula:**  $C_{10}H_{14}N_6O_3$

**Molecular Weight:** 266.25

**Purity:** > 99%

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at +4°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**