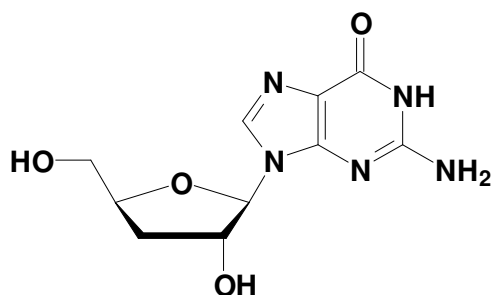


## 3'-Deoxyguanosine

Cat. No.	Amount
N-1048	25 mg



**Cat. No.:** N-1048

**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>4</sub>

**Molecular Weight:** 267.24

**Purity:** > 98%

### 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!**

### Selected References:

Shunji Izuta *et al.* (1996) 3'-Deoxyribonucleotides inhibit Eukaryotic DNA Primase1. *J. Biochem.* **119** (6):1038.

Rasmussen *et al.* (1993) In vivo transcriptional pausing and cap formation on three *Drosophila* heat shock genes. *Proc. Natl. Acad. Sci. USA* **90** 7923

Hamel *et al.* (1976) Interactions of Guanosine Triphosphate Analogues with Elongation Factor G of *Escherichia coli*. *J. Biochem.* **63**:13894.