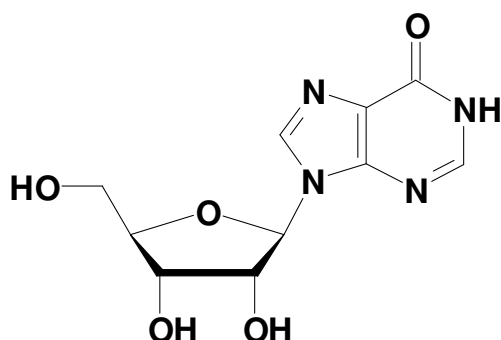


## Inosine

Cat. No.	Amount
N-1049	1 g



**Cat. No.:** N-1049

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>4</sub>O<sub>5</sub>

**Molecular Weight:** 268.22

**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!**