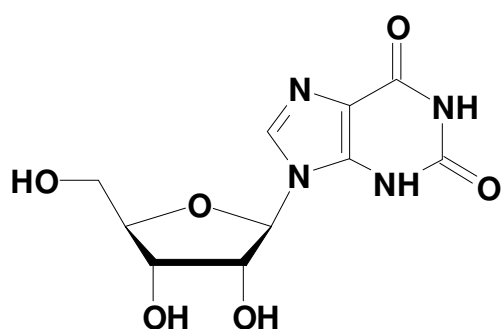


## Xanthosine

Cat. No.	Amount
N-1050	1 g



**Cat. No.:** N-1050

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>4</sub>O<sub>6</sub>

**Molecular Weight:** 284.22

**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!**