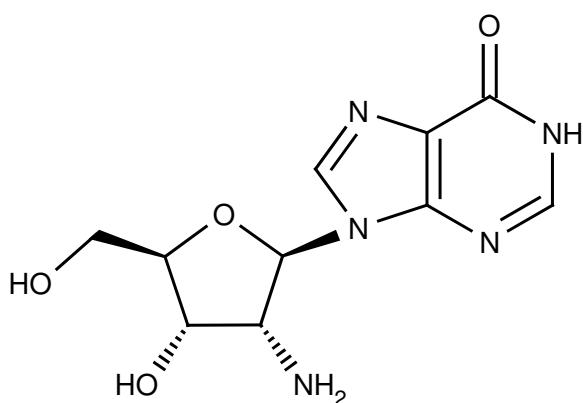


2'-Amino-2'-deoxy-inosine

9-(2-Amino-2-deoxy-β-D-ribofuranosyl)-hypoxanthine

| Cat. No. | Amount |
|--------------|--------|
| N-203-41-250 | 250 mg |
| N-203-41-1G | 1 g |
| N-203-41-10G | 10 g |



Molecular Formula: C₁₀H₁₃N₅O₄

Molecular Weight: 267.24 g/mol

Purity: > 99%

Appearance: white to off-white crystals

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!