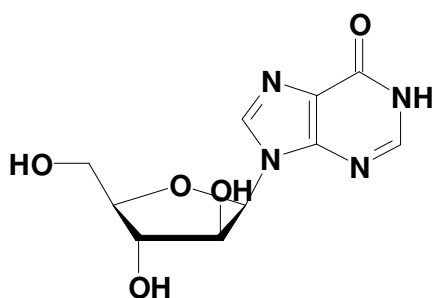


Arabinofuranosyl-Hypoxanthine

9-(β -D-Arabinofuranosyl)-hypoxanthine

Cat. No.	Amount
N-1002	1 g



Cat. No.: N-1002

Molecular Formula: $C_{10}H_{12}N_4O_5$

Molecular Weight: 268.22

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!