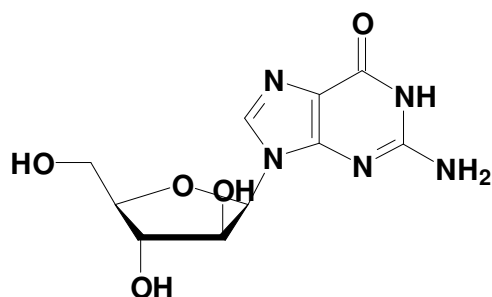


Arabinofuranosyl-Guanine

9-(β -D-Arabinofuranosyl)-guanine

Cat. No.	Amount
N-1003	1 g



Cat. No.: N-1003

Molecular Formula: $\text{C}_{10}\text{H}_{13}\text{N}_5\text{O}_5$

Molecular Weight: 283.24

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!