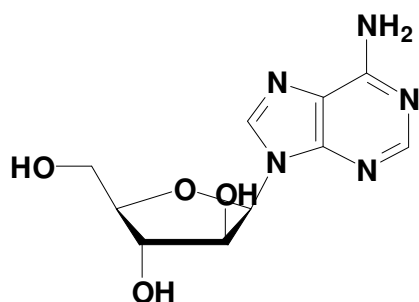


Arabinofuranosyl-Adenine

9-(β -D-Arabinofuranosyl)-adenine

Cat. No.	Amount
N-1005	1 g



Cat. No.: N-1005

Molecular Formula: $C_{10}H_{13}N_5O_4$

Molecular Weight: 267.24

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!