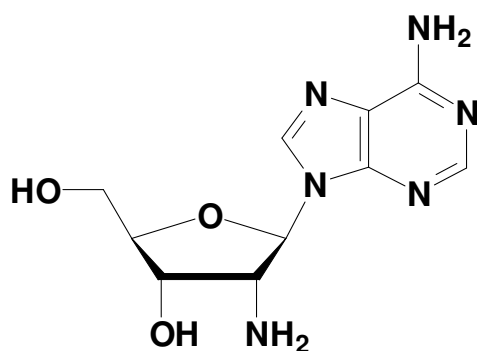


2'-Amino-2'-deoxyadenosine 9-(2-Amino-2-deoxy- β -D-ribofuranosyl)-adenine

Cat. No.	Amount
N-1027	1 g



Cat. No.: N-1027

Molecular Formula: C₁₀H₁₄N₆O₃

Molecular Weight: 266.25

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!