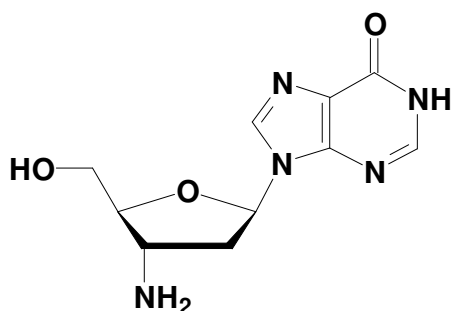


3'-Amino-2',3'-Dideoxyinosine

9-(3-Amino-2,3-dideoxy- β -D-ribofuranosyl)-hypoxanthine

Cat. No.	Amount
N-1009	1 g



Cat. No.: N-1009

Molecular Formula: C₁₀H₁₃N₅O₃

Molecular Weight: 251.24

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!