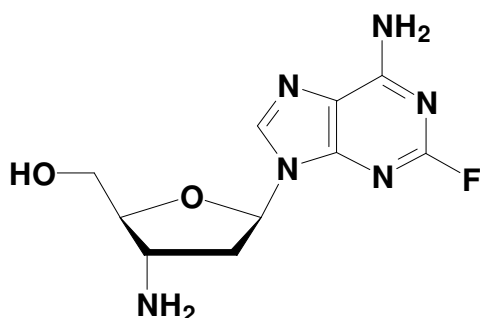


3'-Amino-2',3'-Dideoxy-2-Fluoroadenosine

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

Cat. No.	Amount
N-1301	1 g



Cat. No.: N-1301

Molecular Formula: C₁₀H₁₃FN₆O₂

Molecular Weight: 268.25

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!