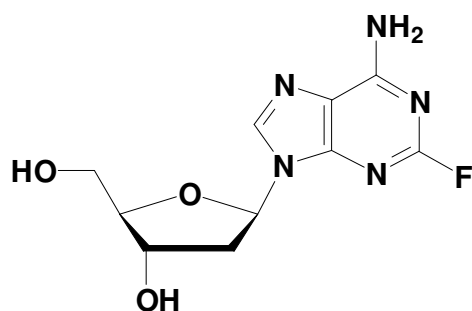


2'-Deoxy-2-fluoroadenosine

9-(2-Deoxy- β -D-ribofuranosyl)-2-fluoroadenine

Cat. No.	Amount
N-1035	1 g



Cat. No.: N-1035

Molecular Formula: C₁₀H₁₂FN₅O₃

Molecular Weight: 269.23

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!