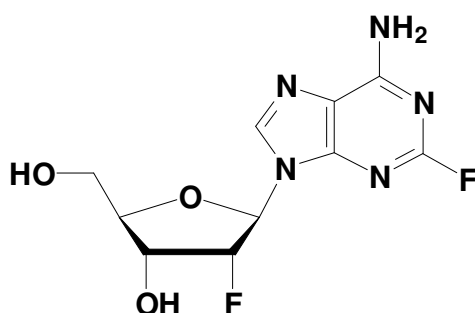


2'-Fluoro-2'-Deoxy-2-fluoroadenosine

9-(2-Fluoro-2-deoxy-β-D-ribofuranosyl)-2-fluoroadenine

Cat. No.	Amount
N-1036	1 g



Cat. No.: N-1036

Molecular Formula: C₁₀H₁₁F₂N₅O₃

Molecular Weight: 287.22

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!