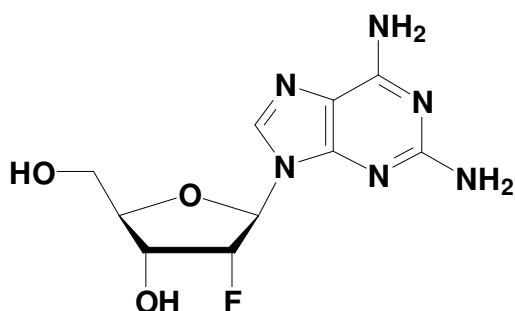


2'-Fluoro-2'-Deoxy-2,6-Diaminopurineriboside

9-(2-Deoxy-2-fluoro-β-D-ribofuranosyl)-2,6-diaminopurine

Cat. No.	Amount
N-1018	1 g



Cat. No.: N-1018

Molecular Formula: C₁₀H₁₃FN₆O₃

Molecular Weight: 284.25

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!