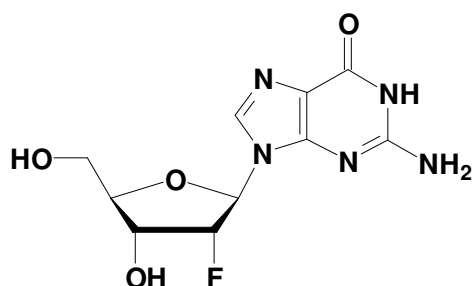


2'-Deoxy-2'-Fluoroguanosine 9-(2-Deoxy-2-fluoro- β -D-ribofuranosyl)-guanine

Cat. No.	Amount
N-1015	1 g



Cat. No.: N-1015

Molecular Formula: C₁₀H₁₂FN₅O₄

Molecular Weight: 285.23

Purity: > 97%

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!