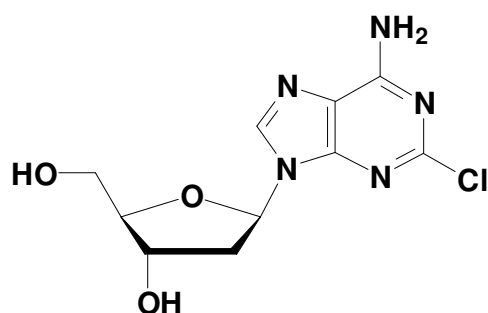


2-Chloro-2'-deoxyadenosine (2-CdA, cladribine)

9-(2-Deoxy- β -D-ribofuranosyl)-2-chloro-adenine

Cat. No.	Amount
N-1025	1 g



Cat. No.: N-1025

Molecular Formula: C₁₀H₁₂ClN₅O₃

Molecular Weight: 285.68

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!