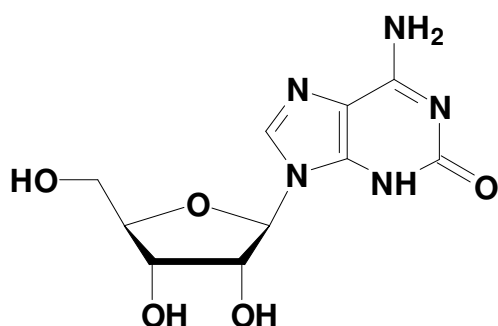


Isoguanosine

9-(β -D-Ribofuranosyl)-isoguanine

Cat. No.	Amount
N-1039	1 g



Cat. No.: N-1039

Molecular Formula: C₁₀H₁₃N₅O₅

Molecular Weight: 283.24

Purity: > 98%

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!