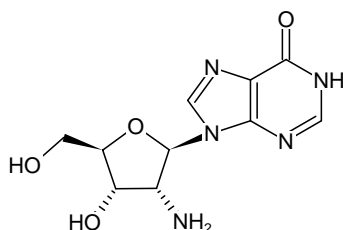


**2'-Amino-2'-deoxyinosine**

9-(2'-Amino-2'-deoxy-β-D-ribofuranosyl)-hypoxanthine

Cat. No.	Amount
N-20341-250	250 mg
N-20341-1G	1 g
N-20341-10G	10 g



Structural formula of 2'-Amino-2'-deoxyinosine

**For research use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at 4 °C**Additional Storage Conditions:** store dark and dry

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 36 months after date of delivery**Molecular Formula:** C<sub>10</sub>H<sub>13</sub>N<sub>5</sub>O<sub>4</sub>**Molecular Weight:** 267.24 g/mol**Purity:** > 99 %**Form:** white to off-white crystals