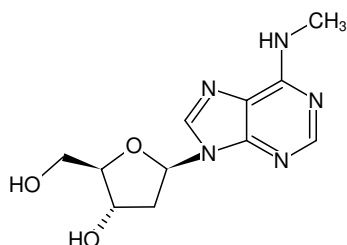


**N<sup>6</sup>-methyl-2'-deoxyadenosine**

Cat. No.	Amount
N-1088	100 mg

Structural formula of N<sup>6</sup>-methyl-2'-deoxyadenosine**For research use only!****Shipping:** shipped at ambient temperature**Storage Conditions:** store at -20 °C**Shelf Life:** 30 months after date of delivery**Molecular Formula:** C<sub>11</sub>H<sub>15</sub>N<sub>5</sub>O<sub>3</sub>**Molecular Weight:** 265.27 g/mol**Exact Mass:** 265.12 g/mol**CAS#:** 2002-35-9**Purity:** ≥ 98 % (HPLC)**Form:** white crystalline powder