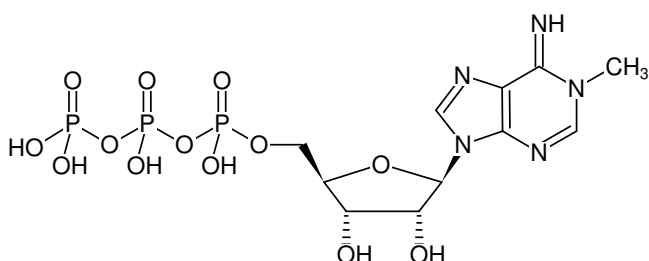


**N<sup>1</sup>-Methyl-ATP - Solid**N<sup>1</sup>-Methyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-1027-1	1 mg
NU-1027-5	5 mg

**Selected References:**Dominissini *et al.* (2016) The dynamic N<sup>1</sup>-methyladenosine methylome in eukaryotic messenger RNA. *Nature* **530 (7591)**:441.Structural formula of N<sup>1</sup>-Methyl-ATP - Solid**For research use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C

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

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>11</sub>H<sub>18</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 521.21 g/mol (free acid)**Exact Mass:** 521.01 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ<sub>max</sub> 257 nm, ε 14.6 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)