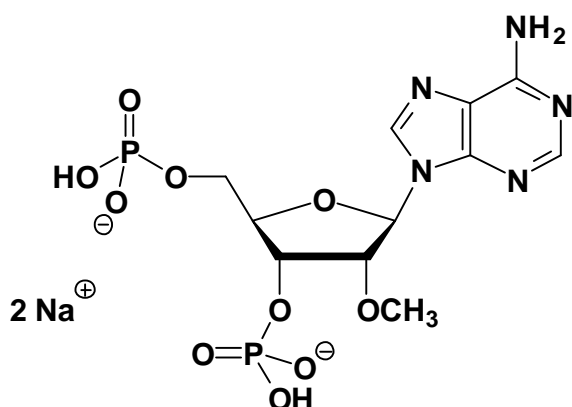


## 2'-OMe-Adenosine-3',5'-bisphosphate (2'-OMe-pAp)

2'-O-Methyl-adenosine-3',5'-bisphosphate, Sodium salt

Cat. No.	Amount
NU-1114S	100 Units
NU-1114L	500 Units



**Cat. No.:** NU-1114

**Molecular Formula:** C<sub>11</sub>H<sub>15</sub>N<sub>5</sub>O<sub>10</sub>P<sub>2</sub> (Anion)

**Molecular Weight:** 439.21 (Anion)

**Purity:** > 95%, clear aqueous solution, pH 7.5

**Storage conditions:**

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**For research use only!**

1 unit = 1 µl of a 10 mM solution