

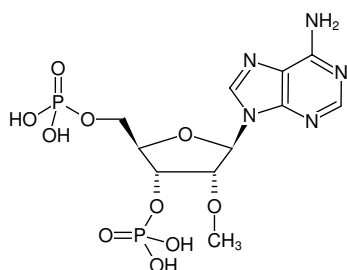


## 2'OMe-Adenosine-3',5'-bisphosphate

2'O-Me-3',5'-pAp

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

Cat. No.	Amount
NU-1114S	100 µl (10 mM)
NU-1114L	5 x 100 µl (10 mM)



Structural formula of 2'OMe-Adenosine-3',5'-bisphosphate

**For general laboratory use.**

**Please centrifuge briefly before opening (volume ≤2 ml).**

**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>17</sub>N<sub>5</sub>O<sub>10</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 441.23 g/mol (free acid)

**Exact Mass:** 441.05 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5

**Spectroscopic Properties:** λ<sub>max</sub> 259 nm, ε 15.3 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)