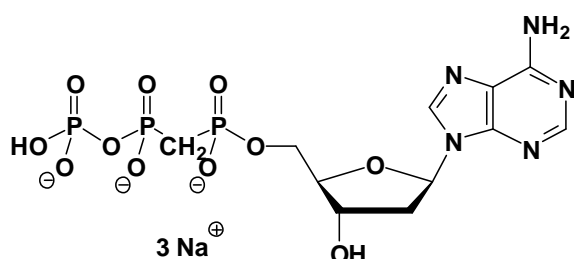


## dApCpp (dAMPCPP)

2'-Deoxy-adenosine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, Sodium salt

Cat. No.	Amount
NU-432S	50 Units
NU-432L	250 Units



**Cat. No.:** NU-432

**Molecular Formula:** C<sub>11</sub>H<sub>15</sub>N<sub>5</sub>O<sub>11</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 486.18 (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  $\mu$ l of a 10 mM solution