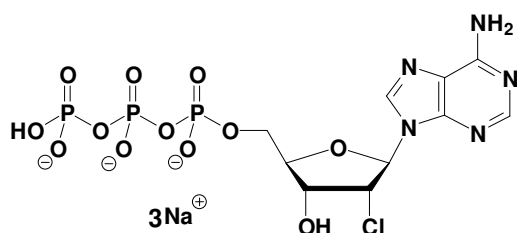


## 2'-Cl-ATP

### 2'-Chloro-Adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-148S	50 Units
NU-148L	250 Units



**Cat. No.:** NU-148

**Molecular Formula:**  $C_{10}H_{12}N_5O_{12}P_3Cl$  (Anion)

**Molecular Weight:** 522.60 (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.

1 unit = 1  $\mu$ l of a 10 mM solution

**For research use only!**