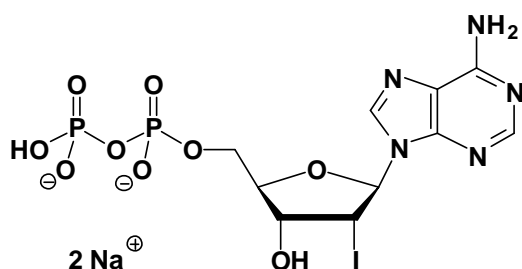


## 2'-I-ADP

2'-Iodo-Adenosine-5'-diphosphate, Sodium salt

Cat. No.	Amount
NU-109S	30 Units
NU-109L	150 Units



**Cat. No.:** NU-109

**Molecular Formula:** C<sub>10</sub>H<sub>12</sub>N<sub>5</sub>O<sub>9</sub>P<sub>2</sub>I (Anion)

**Molecular Weight:** 535.08 (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