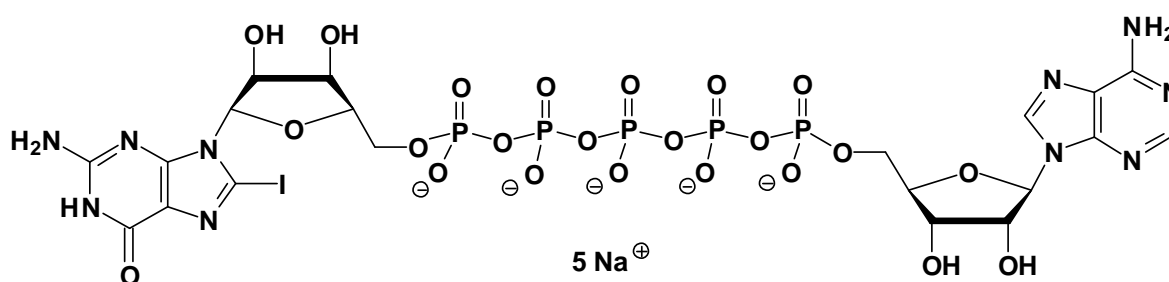


## AP<sub>5</sub>(8I-G)

P<sup>1</sup>-(5'-Adenosyl) P<sup>5</sup>-(5'-(8-Iodo)-guanosyl) pentaphosphate, Sodium salt

Cat. No.	Amount
NU-511S	30 Units
NU-511L	150 Units



**Cat. No.:** NU-511

**Molecular Formula:** C<sub>20</sub>H<sub>23</sub>IN<sub>10</sub>O<sub>23</sub>P<sub>5</sub> (Anion)

**Molecular Weight:** 1052.86 (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