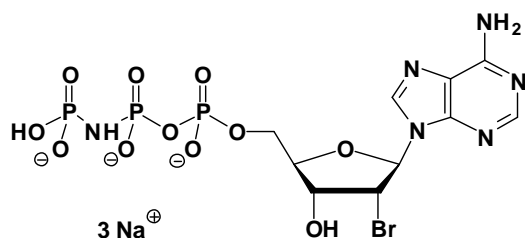


## 2'Br-AppNHp (2'Br-AMPPNP)

2'-Bromo-Adenosine-5'-[ $(\beta,\gamma)$ -imido]triphosphate, Sodium salt

Cat. No.	Amount
NU-104S	20 Units
NU-104L	100 Units



### Selected References:

Gruen *et al.* (1999) 2'-Halo-ATP and -GTP analogues: Rational phasing tools for protein crystallography. *Protein Sci.* **8**:2524.

Gruen *et al.* (1999) Synthesis of 2'-iodo- and 2'-bromo-ATP and GTP analogues as potential phasing tools for X-ray crystallography. *Nucleos. Nucleot.* **18 (1)**:137.

**Cat. No.:** NU-104

**Molecular Formula:** CH<sub>13</sub>N<sub>6</sub>O<sub>11</sub>P<sub>3</sub>Br (Anion)

**Molecular Weight:** 566.07 (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 6 months after date of delivery.

### For research use only!

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