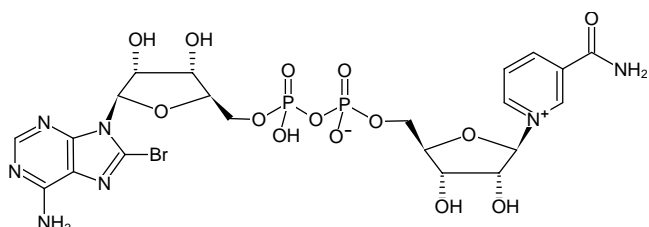


**8-Bromo-AP<sub>2</sub>(Nic)**(8Br-AP<sub>2</sub>(Nic))

Nicotinamide 8-bromo-adenine dinucleotide, Sodium salt

Cat. No.	Amount
NU-523	300 µl (10 mM)

Structural formula of 8-Bromo-AP<sub>2</sub>(Nic)**For research use only!****Shipping:** shipped on blue ice**Storage Conditions:** store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

**Shelf Life:** 12 months after date of delivery**Molecular Formula:** C<sub>21</sub>H<sub>26</sub>BrN<sub>7</sub>O<sub>14</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 742.32 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ±0.5