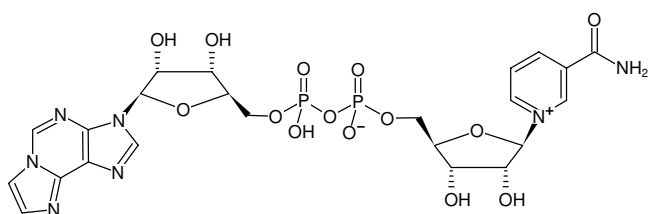



**Etheno-AP<sub>2</sub>(Nic) [ε-AP<sub>2</sub>(Nic)]**

 (Nicotinamide 1,N<sup>6</sup>-ethano-adenine dinucleotide)

 Nicotinamide 1,N<sup>6</sup>-ethano-adenine dinucleotide, Sodium salt

Cat. No.	Amount
NU-522	300 µl (10 mM)


 Structural formula of Etheno-AP<sub>2</sub>(Nic) [ε-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>23</sub>H<sub>27</sub>N<sub>7</sub>O<sub>14</sub>P<sub>2</sub> (free acid)

**Molecular Weight:** 687.45 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solution in water

**Color:** colorless to slightly yellow

**Concentration:** 10 mM - 11 mM

**pH:** 7.5 ±0.5