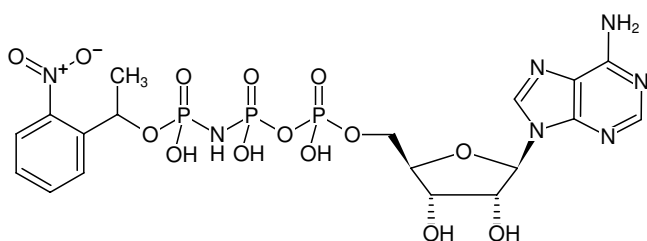


**NPE-caged-AppNHp**

(NPE-caged-AMPPNP)

Adenosine-5'-[( $\beta,\gamma$ )-imido]triphosphate, P<sup>3</sup>-(1-(2-nitrophenyl)-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-305S	20 $\mu$ l (10 mM)
NU-305L	5 x 20 $\mu$ l (10 mM)



Structural formula of NPE-caged-AppNHp

**For research use only!****Shipping:** shipped on gel packs**Storage Conditions:** store at -20 °C**Additional Storage Conditions:** store dark

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

**Shelf Life:** 6 months after date of delivery**Molecular Formula:** C<sub>18</sub>H<sub>24</sub>N<sub>7</sub>O<sub>14</sub>P<sub>3</sub> (free acid)**Molecular Weight:** 655.35 g/mol (free acid)**Exact Mass:** 655.06 g/mol (free acid)**CAS#:** 116271-21-7**Purity:**  $\geq$  95 % (HPLC)**Form:** solution in water**Color:** colorless to slightly yellow**Concentration:** 10 mM - 11 mM**pH:** 7.5  $\pm$  0.5**Spectroscopic Properties:**  $\lambda_{\max}$  260 nm,  $\epsilon$  18.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)**Applications:**

Agonistic ligand, mainly for nucleoside receptor A<sub>1</sub>. Nucleosidephosphates stabilized against hydrolytic degradation can directly bind to nucleoside receptors. The caged form is protected during uptake and transport in animal experiments and can be well-directed released through activation at the target tissue.

**Selected References:**

Volonte *et al.* (2009) Membrane components and purinergic signalling: the purinome, a complex interplay among ligands, degrading enzymes, receptors and transporters. *FEBS J.* **276**:318.

Yegutkin (2008) Nucleotide and nucleoside converting enzymes: Important modulators of purinergic signalling cascade. *Biochim. Biophys. Acta* **1783**:673.

Williams *et al.* (1986) Effects of purine nucleotides on the binding of [<sup>3</sup>H]cyclopentyladenosine to adenosine A<sub>1</sub>-receptors in rat brain membranes. *J. Neurochem.* **47** (1):88.