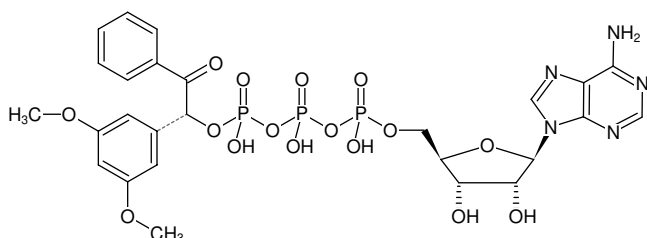


**DMB-caged-ATP**Adenosine-5'-triphosphate, P³-(1-(3',5'-dimethoxyphenyl)-2-oxo-2-phenyl-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-309S	10 µl (10 mM)
NU-309L	5 x 10 µl (10 mM)



Structural formula of DMB-caged-ATP

For research use only!**Shipping:** shipped on blue ice**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: 12 months after date of delivery**Molecular Formula:** C₂₆H₃₀N₅O₁₆P₃ (free acid)**Molecular Weight:** 761.46 g/mol (free acid)**Exact Mass:** 761.09 g/mol (free acid)**CAS#:** 159899-51-1**Purity:** ≥ 95 % (HPLC)**Form:** clear aqueous solution**Concentration:** 10 mM - 11 mM**pH:** 7.5 ± 0.5**Spectroscopic Properties:** λ_{max} 256 nm, ε 25.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)**Applications:****Ligand for purinergic receptors:**

The nucleotide can be transported extra- or intracellular in a protected form to the target. After activation by well-defined conditions the liberated ATP can interact with P2X- and P2Y-receptors. Interacting subreceptor types and corresponding references are listed in Data sheet #NU-1010.

Agonistic ligand, mainly for nucleoside receptor A₁

Nucleoside-triphosphates can be converted by different membrane-bound phosphatases into nucleosides acting as nucleoside receptor ligands. 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.

Apell *et al.* (1998) Partial reactions of the Na,K-ATPase: kinetic analysis and transport properties. *Acta Physiol. Scand.* **163**:235.

Sokolov *et al.* (1998) Fast transient currents in Na,K-ATPase induced by ATP concentration jumps from the P-3-[1-(3',5'-dimethoxyphenyl)-2-oxo]ethyl ester of ATP. *Biophys. J.* **74** (5):2285.

Sokolov *et al.* Fast transient currents in the Na,K-ATPase induced by ATP concentration jump experiments from DMBcaged ATP. *Biophys. J.* **72** (2):242.

Thirlwell *et al.* (1995) Inhibition of unloaded shortening velocity in permeabilized muscle-fibers by caged-ATP compounds. *J. Muscle Res. Cell M.* **16** (2):131.

Thirlwell *et al.* (1994) Kinetics of relaxation from rigor of permeabilized fast-twitch skeletal fibers from the rabbit using a novel caged-ATP and apyrase. *Biophys. J.* **67** (6):2436.

Corrie *et al.* (1992) Synthetic, mechanistic and photochemical studies of phosphate-esters of substituted benzoines. *J. Chem. Soc. Perkin Trans.* **1** (18):2409.

Corrie *et al.* (1992) The development and application of photosensitive caged compounds to aid time-resolved structure determination of macromolecules. *Philos. T. Roy. Soc. A* **340** (1657):233.