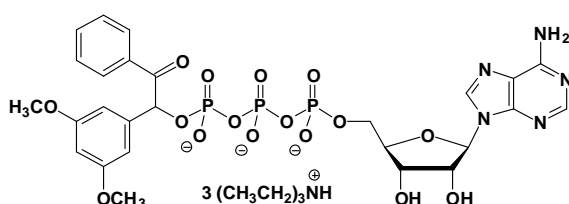


DMB-caged-ATP

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

Cat. No.	Amount
NU-309S	10 Units
NU-309L	50 Units



Cat. No.: NU-309

Molecular Formula: C₂₆H₂₇N₅O₁₆P₃ (Anion)

Molecular Weight: 758.44 (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 12 months after date of delivery. Store in a dark place and protect from light!

For research use only!

1 unit = 1 µl of a 10 mM solution

Selected References:

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-phenyl-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 benzoin. *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