

**BzBzATP (BzATP)**

2'/3'-O-(4-Benzoylbenzoyl)adenosine-5'-triphosphate, Tri(triethylammonium) salt

Cat. No.	Amount
NU-1620-5	5 mg
NU-1620-25	25 mg

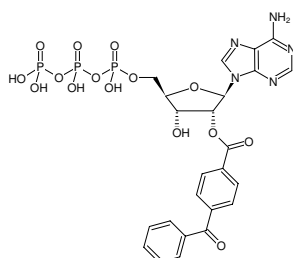
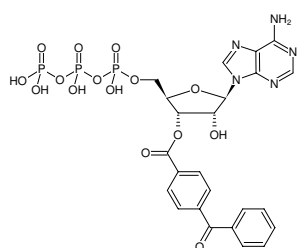
Specific Ligands:Agonistic ligand for P2X₇ receptor^[1,2]**Note : Product is hygroscopic!****Selected References:**

[1] Gibbons *et al.* (2001) P2X₇ receptors in rat parotid acinar cells: formation of large pores. *J. Auton. Pharmacol.* **21** (4):181.

[2] Huong-Zhen Hu *et al.* (2001) P2X₇ receptors in the enteric nervous system of guinea pig small intestine. *J. Comp. Neurol.* **440** (3):299.

Miras-Portugal *et al.* (2007) P2X agonist BzATP interferes with amplex-red-coupled fluorescence assays. *Anal. Biochem.* **367**:140.

Mahmood *et al.* (1989) Serine-324 of myosin's heavy chain is photoaffinity-labeled by 3' (2')-O- (4-benzoylbenzoyl)adenosine triphosphate. *Biochemistry* **28**:3989.



Structural formula of BzBzATP (BzATP)

For general laboratory use.**Shipping:** shipped on gel packs**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:** C₂₄H₂₄N₅O₁₅P₃*3(C₆H₁₅N) [tri(triethylammonium salt)]; C₂₄H₂₄N₅O₁₅P₃ (free acid)**Molecular Weight:** 1018.97 g/mol [tri(triethylammonium salt)]; 715.40 g/mol (free acid)**Exact Mass:** 715.05 g/mol (free acid)**CAS#:** 112898-15-4**Purity:** ≥ 95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:** λ_{max} 261 nm, ε 41.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)