

**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

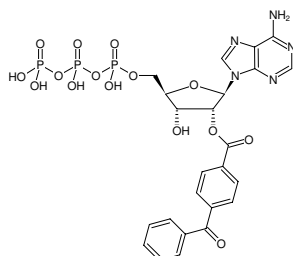
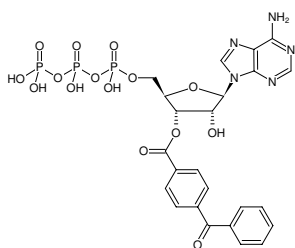
**Specific Ligands:**Agonistic ligand for P2X<sub>7</sub> receptor<sup>[1,2]</sup>**Note : Product is hygroscopic!****Selected References:**

[1] Gibbons *et al.* (2001) P2X<sub>7</sub> receptors in rat parotid acinar cells: formation of large pores. *J. Auton. Pharmacol.* **21** (4):181.

[2] Huang-Zhen Hu *et al.* (2001) P2X<sub>7</sub> 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 research use only!****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:** C<sub>24</sub>H<sub>24</sub>N<sub>5</sub>O<sub>15</sub>P<sub>3</sub>\*3(C<sub>6</sub>H<sub>15</sub>N) [tri(triethylammonium salt)]; C<sub>24</sub>H<sub>24</sub>N<sub>5</sub>O<sub>15</sub>P<sub>3</sub> (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<sub>max</sub> 261 nm, ε 41.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)