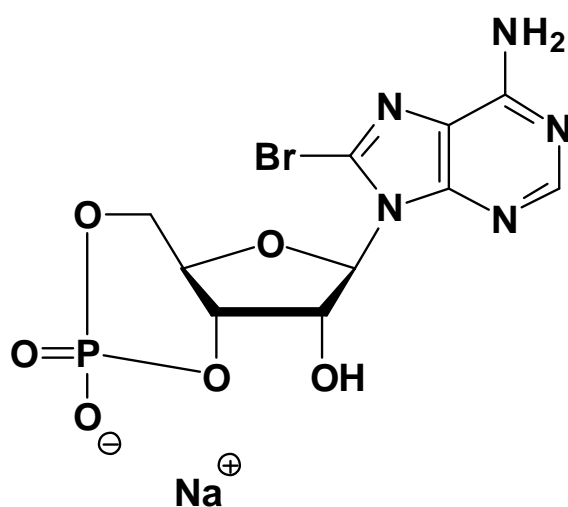


## 8Br-cAMP

8-Bromo-Adenosine-3',5'-cyclic monophosphate, Sodium salt

Cat. No.	Amount
NU-1502S	10 mg
NU-1502L	50 mg



**Cat. No.:** NU-1502

**Molecular Formula:** C<sub>10</sub>H<sub>10</sub>N<sub>5</sub>O<sub>6</sub>PBr (Anion)

**Molecular Weight:** 407.09 (Anion)

**Purity:** > 95%, lyophilized solid

### 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.

**For research use only!**

### Selected References:

Paz *et al.* (1999) Corticotropin increases protein tyrosine phosphatase activity by a cAMP-dependent mechanism in rat adrenal gland. *Eur. J. Biochem.* **265**:911.

Martin *et al.* (1999) Comparison of expression and regulation of the high-density lipoprotein receptor SR-BI and the lowdensity lipoprotein receptor in human adrenocortical carcinoma NCI-H295 cells. *Eur. J. Biochem.* **261**:481.

Chen *et al.* (1998) Up-regulation of the cAMP/PKA pathway inhibits proliferation, induces differentiation, and leads to apoptosis in malignant gliomas. *Lab. Invest.* **78**:165.

Hei *et al.* (1991) Lack of correlation between activation of cyclic AMP-dependent protein kinase and inhibition of contraction of rat vas deferens by cyclic AMP analogs. *Mol. Pharmacol.* **39**:233.

Malgaroli *et al.* (1989) Control of cytosolic free calcium in rat and chicken osteoclasts. The role of extracellular calcium and calcitonin. *J. Biol. Chem.* **264**:14342.

Schubart *et al.* (1977) Cyclic adenosine 3':5'-monophosphate-mediated insulin secretion and ribosomal protein phosphorylation in a hamster islet cell tumor. *J. Biol. Chem.* **252**:92.

Martin *et al.* (1981) Growth inhibition by adenosine 3',5'-monophosphate derivatives does not require 3',5' phosphodiester linkage. *Science* **213**:1120.

Meyer *et al.* (1974) Analogs of cyclic AMP and cyclic GMP: general methods of synthesis and the relationship of structure to enzymic activity. *Life Sci.* **14**:1019.