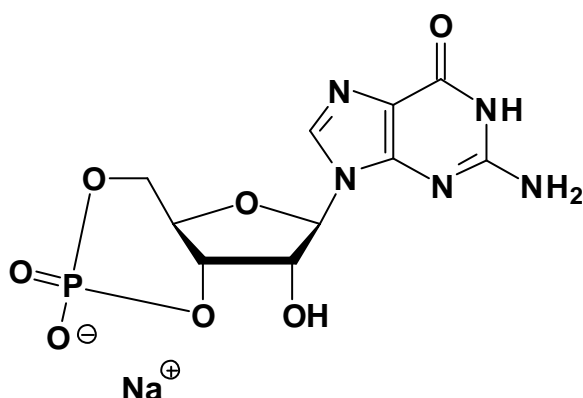


cGMP

Guanosine-3',5'-cyclic monophosphate, Sodium salt

Cat. No.	Amount
NU-1501S	50 mg
NU-1501L	250 mg



Cat. No.: NU-1501

Molecular Formula: C₁₀H₁₁N₅O₇P (Anion)

Molecular Weight: 344.2 (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:

Nisoli et al. (2003) Mitochondrial biogenesis in mammals: the role of endogenous nitric oxide. *Science* **299** (5608):896.

Ahumada et al. (2003) Signaling of rat Frizzled-2 through phosphodiesterase and cyclic GMP. *Science* **298** (5600):2006.

Rybalkin et al. (2003) PDE5 is converted to an activated state upon cGMP binding to the GAF A domain. *EMBO J.* **22** (3):469.

Butt et al. (2003) Actin binding of human LIM and SH3 protein (LASP) is regulated by cGMP- and cAMP-dependent protein kinase phosphorylation on Ser-146. *J. Biol. Chem.* **2003 Feb 5** [epub ahead of print].

Leckie et al. (2003) The NO pathway acts late during the fertilization response in sea urchin eggs. *J. Biol. Chem.* **2003 Jan 22** [epub ahead of print]

Ludidi et al. (2003) Identification of a Novel Protein with Guanylyl Cyclase Activity in *Arabidopsis thaliana*. *J. Biol. Chem.* **278** (8):6490.

Coates et al. (2002) Antagonistic pathways in neurons exposed to body fluid regulate social feeding in *Caenorhabditis elegans*. *Nature* **419** (6910):925.

Bosgraaf et al. (2002) A novel cGMP signalling pathway mediating myosin phosphorylation and chemotaxis in *Dictyostelium*. *EMBO J.* **21** (17):4560.

Higgins et al. (2002) Molecular architecture of a retinal cGMP-gated channel: the arrangement of the cytoplasmic domains. *EMBO J.* **21** (9):2087.

Castellani et al. (2002) Cis and trans interactions of L1 with neuropilin-1 control axonal responses to semaphorin 3A. *EMBO J.* **21** (23):6348.

Park et al. (2002) Nitric Oxide-Cyclic GMP Signaling Pathway in the Regulation of Rabbit Clitoral Caverosum Tone. *Exp. Biol. Med.* (Maywood) **227** (11):1022.