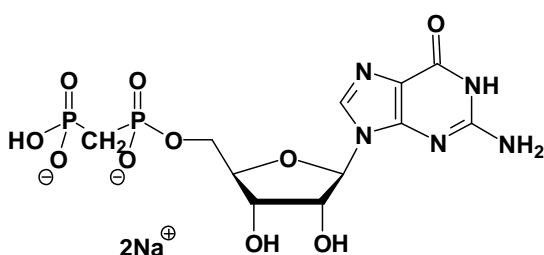


GpCp (GMPCP)

Guanosine-5'-[(α,β)-methyleno]diphosphate, Sodium salt

Cat. No.	Amount
NU-414S	150 Units
NU-414L	750 Units



Cat. No.: NU-414

Molecular Formula: C₁₁H₁₅N₅O₁₀P₂ (Anion)

Molecular Weight: 439.21 (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.

For research use only!

1 unit = 1 μ l of a 10 mM solution

Selected References:

Muller-Reichert *et al.* (1998) Structural changes at microtubule ends accompanying GTP hydrolysis: information from a slowly hydrolyzable analogue of GTP, guanylyl (α,β)methylenediphosphonate. *Proc. Natl. Acad. Sci. USA* **95** (7):3661.

Vulevic *et al.* (1997) Role of guanine nucleotides in the vinblastine-induced self-association of tubulin: effects of guanosine α,β -methylenetriphosphate and guanosine α,β -methylenediphosphate. *Biochemistry-US* **36** (42):12828.

Hyman *et al.* (1995) Structural-changes accompanying GTP hydrolysis in microtubules - information from a slowly hydrolyzable analog guanylyl-(α,β)-methylene-diphosphonate. *J. Cell Biol.* **128** (1-2):117.

Hyman *et al.* (1992) Role of GTP hydrolysis in microtubule dynamics - information from a slowly hydrolyzable analog, GMPCPP. *Mol. Biol. Cell* **3** (10):1155.