



pGpG

P¹-(5'-Guanosyl)-P²-(5'-guanosyl)-(3'→5')-diphosphate, Sodium salt





Structural formula of pGpG

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: C₂₀H₂₆N₁₀O₁₅P₂ (free acid)

Molecular Weight: 708.43 g/mol (free acid)

Exact Mass: 708.11 g/mol (free acid)

CAS#: 33008-99-0

Purity: ≥ 95 % (HPLC)

Form: solid

Color: white to off-white

Spectroscopic Properties: λ_{max} 252 nm, ϵ 25.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Description:

The linear diguanylate pGpG is the first phosphodiesterase-catalyzed hydrolysis product of cylic di-GMP (c-diGMP). C-diGMP has been identified as a novel second messenger in bacteria that triggers various physiological changes such as virulence gene expression or cell differentiation^[1-5]. The role of pGpG in c-diGMP signalling is under investigation.

Selected References:

[1] Roemling et al. (2005) C-diGMP: the Dawning of a novel bacterial singalling system. Mol. Microbiol. 57:629.

[2] Hengge (2009) Principles of c-di-GMP signalling in bacteria. Nat. Rev. Microbiol. 7 (4):263.

[3] Sudarsan et al. (2008) Riboswitches in Eubacteria Sense the Second Messenger cyclic di-GMP. Science 321:411.

[4] Rao et al. (2010) YybT is a Signaling Protein that Contains a Cyclic Dinucleotide Phosphodiesterase Domain and a GGDEF Domain with ATPase Activity. J. Biol. Chem. 285:473.

[5] Christen et al. (2007) DgrA is a Member of a New Family of Cyclic Diguanosine Monophosphate Receptors and Controls Flagellar Motor Function in Caulobacter crescentus. Proc. Nat. Acad. Sci. USA 104:4112.

