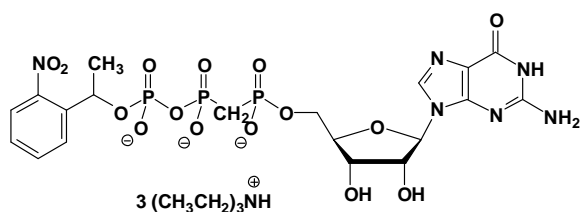


## NPE-caged-GpCpp (NPE-caged-GMPCPP)

Guanosine-5'-[( $\alpha,\beta$ )-methylene]triphosphate, P<sup>3</sup>-(1-(2-nitrophenyl)-ethyl)-ester, Triethylammonium salt

Cat. No.	Amount
NU-306S	10 Units
NU-306L	50 Units



**Cat. No.:** NU-306

**Molecular Formula:** C<sub>19</sub>H<sub>22</sub>N<sub>6</sub>O<sub>15</sub>P<sub>3</sub> (Anion)

**Molecular Weight:** 667.33 (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. Store in a dark place and protect from light!

### For research use only!

1 unit = 1  $\mu$ l of a 10 mM solution