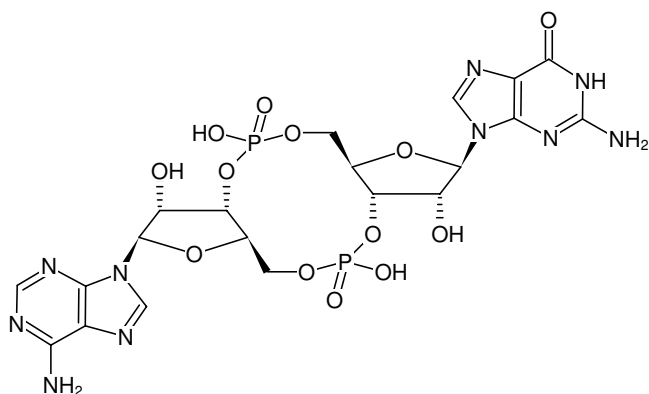


**3',3'-cGAMP**

3',3'-cyclic GMP-AMP

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

Cat. No.	Amount
NU-986S	1 $\mu$ mol
NU-986L	5 x 1 $\mu$ mol



Structural formula of 3',3'-cGAMP

**For research use only!****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:** C<sub>20</sub>H<sub>24</sub>N<sub>10</sub>O<sub>13</sub>P<sub>2</sub> (free acid)**Molecular Weight:** 674.41 g/mol (free acid)**Exact Mass:** 674.10 g/mol (free acid)**Purity:**  $\geq$  95 % (HPLC)**Form:** solid**Color:** white to off-white**Spectroscopic Properties:**  $\lambda_{\max}$  260 nm,  $\epsilon$  25.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)