

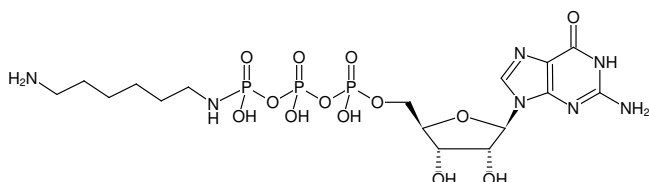


## $\gamma$ -[(6-Aminoethyl)-imido]-GTP

$\gamma$ -[(6-Aminoethyl)-imido]-guanosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-816-1	1 mg
NU-816-5	5 mg

**Please note:** This compound contains a phosphoramidate linkage which is hydrolyzed at pH <7.0.



Structural formula of  $\gamma$ -[(6-Aminoethyl)-imido]-GTP

### 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:** C<sub>16</sub>H<sub>30</sub>N<sub>7</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 621.37 g/mol (free acid)

**Exact Mass:** 621.11 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:**  $\lambda_{\text{max}}$  252 nm,  $\epsilon$  13.7 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)