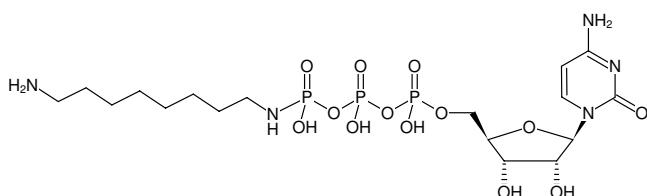


 **$\gamma$ -[(8-Aminoethyl)-imido]-CTP** $\gamma$ -[(8-Aminoethyl)-imido]-cytidine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-844-5	5 mg
NU-844-25	25 mg

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



Structural formula of  $\gamma$ -[(8-Aminoethyl)-imido]-CTP

**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>17</sub>H<sub>34</sub>N<sub>5</sub>O<sub>13</sub>P<sub>3</sub> (free acid)

**Molecular Weight:** 609.40 g/mol (free acid)

**Exact Mass:** 609.14 g/mol (free acid)

**Purity:** ≥ 95 % (HPLC)

**Form:** solid

**Color:** white to off-white

**Spectroscopic Properties:**  $\lambda_{\max}$  271 nm,  $\epsilon$  8.9 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)