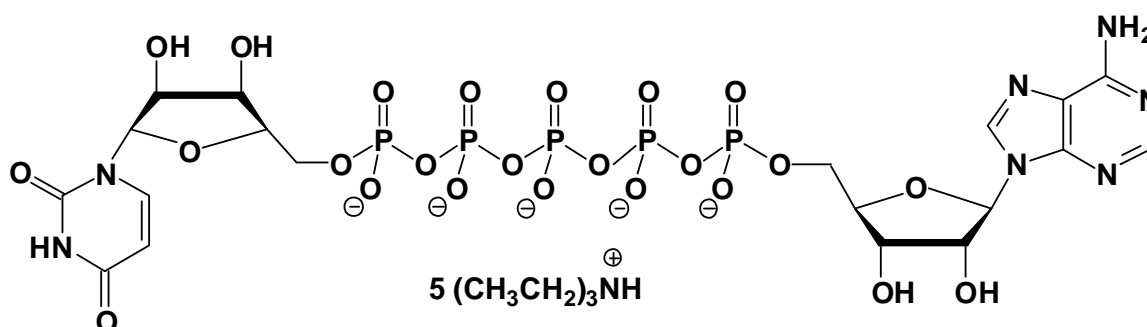


## AP<sub>5</sub>U

P<sup>1</sup>-(5'-Adenosyl) P<sup>5</sup>-(5'-uridyl) pentaphosphate, Sodium salt

Cat. No.	Amount
NU-505S	50 Units
NU-505L	250 Units



**Cat. No.:** NU-505

**Molecular Formula:** C<sub>19</sub>H<sub>23</sub>N<sub>7</sub>O<sub>24</sub>P<sub>5</sub> (Anion)

**Molecular Weight:** 888.29 (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.

**For research use only!**

1 unit = 1 μl of a 10 mM solution

**Selected References:**

Cheng *et al.* (1986) Homogeneous uridine kinase from Ehrlich ascites tumor: substrate specificity and inhibition by bisubstrate analogs. *Mol Pharmacol.* **30** (2):159.