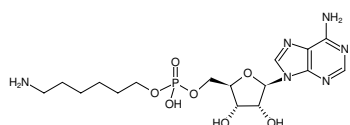


 **$\alpha$ -(6-Aminoheptyl)-AMP** $\alpha$ -(6-Aminoheptyl)-adenosine-5'-monophosphate, Sodium salt

Cat. No.	Amount
NU-859-1	1 mg
NU-859-5	5 mg

Structural formula of  $\alpha$ -(6-Aminoheptyl)-AMP**For research use only!****Shipping:** shipped on blue ice**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>27</sub>N<sub>6</sub>O<sub>7</sub>P (free acid)**Molecular Weight:** 446.40 g/mol (free acid)**Exact Mass:** 446.17 g/mol (free acid)**Purity:** ≥ 95 % (HPLC)**Form:** solid