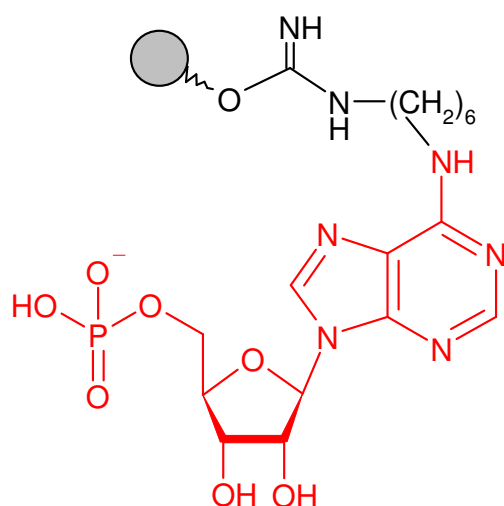


## N<sup>6</sup>-(6-Amino)hexyl-AMP - Sepharose<sup>®</sup>

Cat. No.	Amount
AC-145S	1 ml (bulk material)
AC-145L	5 ml (bulk material)
AC-145C	1 ml (syringe column)
AC-145SC	0.2 ml (scr. column)
AC-145MP	1 ml (MPLC column)



N<sup>6</sup>-(6-Amino)hexyl-adenosine-5'-monophosphate, N<sup>6</sup>-(6-Amino)hexyl-AMP immobilized on Agarose

**Degree of substitution:** 5 μmol N<sup>6</sup>-(6Amino)-hexyl-AMP/ml Gel

**Linker:** none

**Storage buffer:** 20% Ethanol

**Storage:** Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Long term storage at 4°C.

**Expiry:** 12 month

® Sepharose is a trademark of Amersham Biosciences.