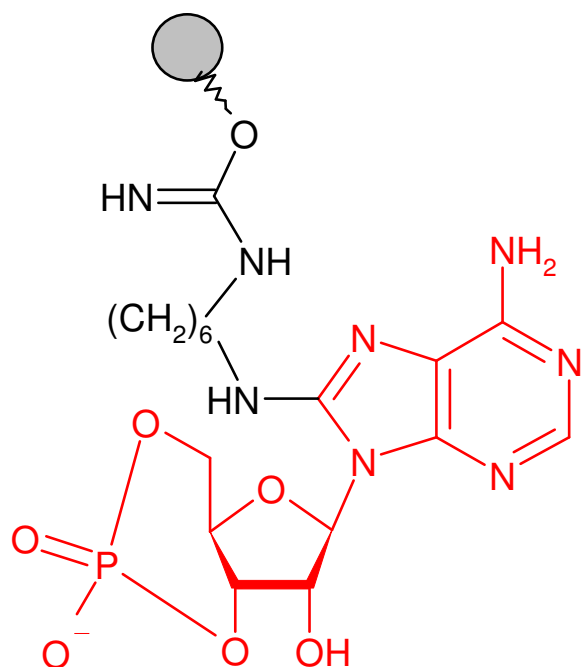


8-Amino-hexyl-cAMP - Sepharose®

Cat. No.	Amount
AC-146S	1 ml (bulk material)
AC-146L	5 ml (bulk material)
AC-146C	1 ml (syringe column)
AC-146SC	0.2 ml (scr. column)
AC-146MP	1 ml (MPLC column)



(8-[(6-Amino)hexyl]amino-adenosine-3', 5'-cyclic-mono-phosphate, 8-[(6-Amino)hexyl]amino-cAMP immobilized on Agarose

Degree of substitution: $5 \mu\text{mol}$ 8-[(6-Amino)hexyl]amino-cAMP/ml Gel $5 \mu\text{mol}$ 2'/3'-EDA-m7-GDP/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.