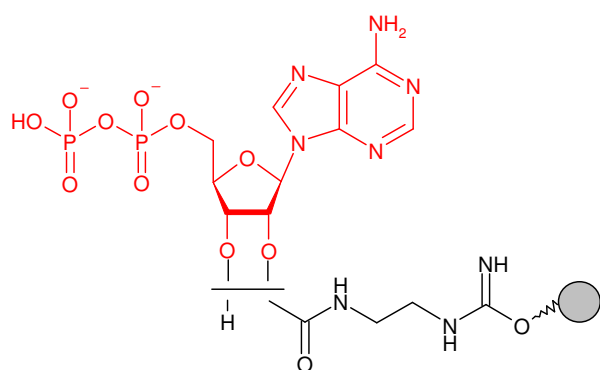


2'/3'-EDA-ADP - Sepharose®

Cat. No.	Amount
AC-144S	1 ml (bulk material)
AC-144L	5 ml (bulk material)
AC-144C	1 ml (syringe column)
AC-144SC	0.2 ml (scr. column)
AC-144MP	1 ml (MPLC column)



2'/3'-O-(2-Aminoethyl-carbamoyl)-Adenosine -5'-
diphosphate (2'/3'-EDA-ADP) immobilized on Agarose

Degree of substitution: 5 μmol 2'/3'-EDA-ADP/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.