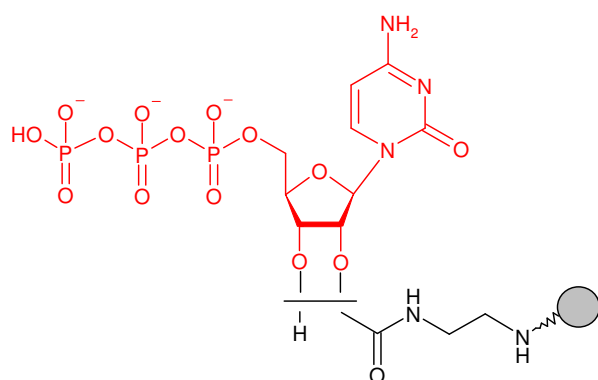


## 2'/3'-EDA-CTP - Sepharose®

Cat. No.	Amount
AC-134S	1 ml (bulk material)
AC-134L	5 ml (bulk material)
AC-134C	1 ml (syringe column)
AC-134SC	0.2 ml (scr. column)
AC-134MP	1 ml (MPLC column)



2'/3'-EDA-Cytidine 5'-triphosphate (2'/3'-EDA-CTP, 2'/3'-O-(2-Aminoethyl-carbamoyl)-cytidine 5'-triphosphate) immobilized on Agarose, suitable for purification of CTP-binding proteins.

**Degree of substitution:** 5  $\mu$ mol 2'/3'-EDA-CTP/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.