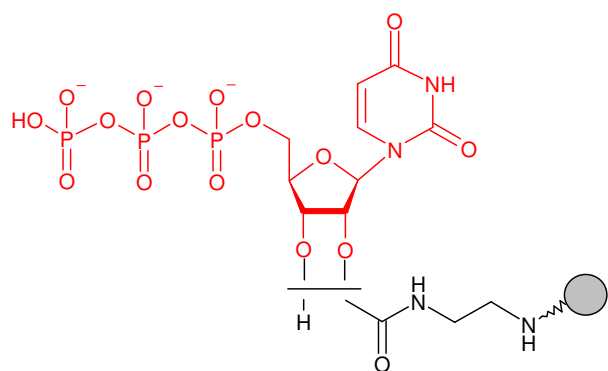


2'/3'-EDA-UTP - Sepharose®

Cat. No.	Amount
AC-133S	1 ml (bulk material)
AC-133L	5 ml (bulk material)
AC-133C	1 ml (syringe column)
AC-133SC	0.2 ml (scr. column)
AC-133MP	1 ml (MPLC column)



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

Degree of substitution: 5 μ mol 2'/3'-EDA-UTP/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.