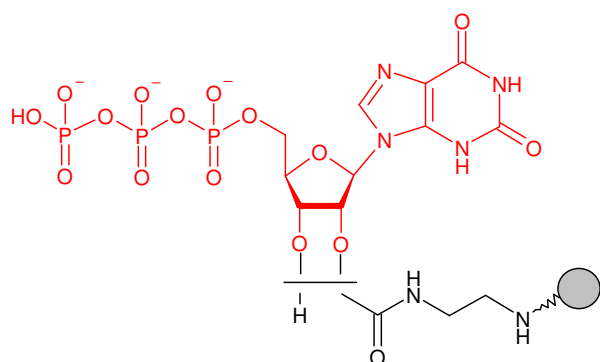


2'/3'-EDA-XTP - Sepharose®

Cat. No.	Amount
AC-136S	1 ml (bulk material)
AC-136L	5 ml (bulk material)
AC-136C	1 ml (syringe column)
AC-136SC	0.2 ml (scr. column)
AC-136MP	1 ml (MPLC column)



2'/3'-EDA-xanthosine 5'-triphosphate (2'/3'-EDA-XTP, 2'/3'-O-(2-Aminoethyl-carbamoyl)-Xan-thosine 5'-triphosphate) immobilized on Agarose, suitable for purification of XTP-binding pro-teins.

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