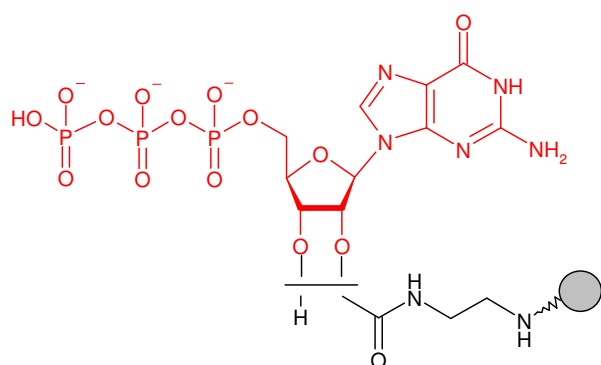


## 2'/3'-EDA-GTP - Sepharose®

Cat. No.	Amount
AC-132S	1 ml (bulk material)
AC-132L	5 ml (bulk material)
AC-132C	1 ml (syringe column)
AC-132SC	0.2 ml (scr. column)
AC-132MP	1 ml (MPLC column)



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

**Degree of substitution:** 5  $\mu\text{mol}$  2'/3'-EDA-GTP/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.