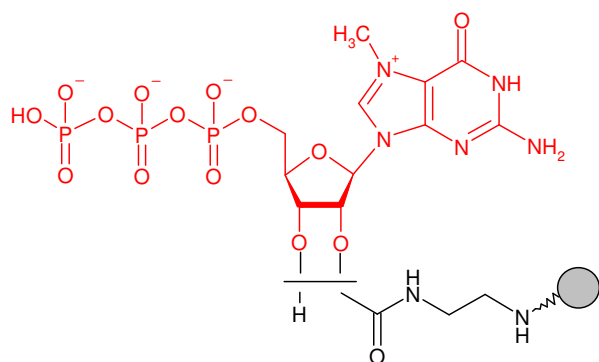


2'/3'-EDA-m⁷GTP - Sepharose[®]

Cat. No.	Amount
AC-142S	1 ml (bulk material)
AC-142L	5 ml (bulk material)
AC-142C	1 ml (syringe column)
AC-142SC	0.2 ml (scr. column)
AC-142MP	1 ml (MPLC column)



2'/3'-EDA-7-Methylguanosine-5'-triphosphate (2'/3'-EDA-m⁷GTP) immobilized on Agarose, suitable for purification of eukaryotic mRNA Cap-Binding Protein.

Degree of substitution: 5 μ mol 2'/3'-EDA-m⁷-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.

Selected References:

Webb *et al.* (1984) Purification of the mes-senger RNA cap-binding protein using a new affinity medium *Biochemistry*. **17;23(2)**:177-81.

Morino *et al.* (1996) Crystallization and pre-liminary X-ray diffraction study of recombi-nant human eukaryotic initiation factor-4E. *J Biochem (Tokyo)*. **119(2)**:224-5.