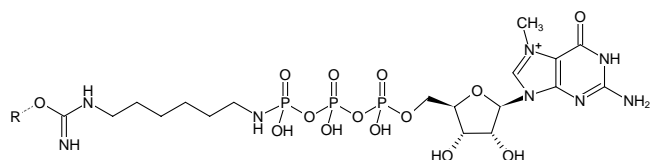




Immobilized γ -Aminoethyl- m^7 GTP

7-Methyl-guanosine 5'-triphosphate (m^7 GTP) immobilized on Agarose
 γ -Aminoethyl- m^7 GTP-Agarose

Cat. No.	Amount
AC-141S	1 ml
AC-141L	5 ml



Structural formula of Immobilized γ -Aminoethyl- m^7 GTP

	Agarose characteristics
Bead/Particle size	45-165 μ m
Recommended linear flow rate	11.5 cm/h
Maximum pressure	0.25 bar (3.6 psi)
pH stability	short term: 4 - 9 / long term: 7.5
Chemical stability	Stable to all solutions commonly used in gel filtration including 8 M urea and 6 M guanidine hydrochloride Not stable in organic solvents!
Sterilization	Not autoclavable!

R= Agarose

For research use only!

Shipping: shipped at 4 °C

Storage Conditions: store at 4 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Degree of substitution: 5 μ mol - 7 μ mol m^7 GTP/ml gel

Storage buffer: 20% Ethanol

Please note:

For the purification of eukaryotic mRNA cap-binding proteins we recommend AC-155. AC-141 contains a phosphoramidate linkage which is hydrolyzed at pH < 7.

Selected References:

Marti *et al.* (2017) No Escaping the Rat Race: Simulated Night Shift Work Alters the Time-of-Day Variation in BMAL1 Translational Activity in the Prefrontal Cortex. *Front Neural Circuits*. doi:10.3389.