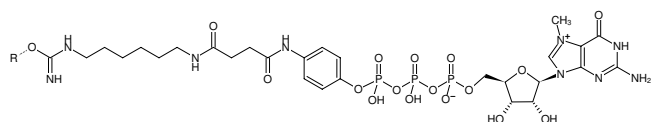




Immobilized γ -Aminophenyl- m^7 GTP (C_{10} -spacer)

γ -Aminophenyl- m^7 GTP (C_{10} -spacer)-Agarose

Cat. No.	Amount
AC-155S	1 ml
AC-155L	5 ml



Structural formula of Immobilized γ -Aminophenyl- m^7 GTP (C_{10} -spacer)

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 general laboratory use.

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

Applications:

Suitable for purification of eukaryotic mRNA cap-binding proteins.

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

Storage buffer: 20% Ethanol, 0.1 M Tris-HCl pH 7.5

Selected References:

Coots *et al.* (2017) m6A Facilitates eIF4F-Independent mRNA Translation. *Mol. Cell.* doi:10.1016.

Hou *et al.* (2017) PERK Signal-Modulated Protein Translation Promotes the Survivability of Dengue 2 Virus-Infected Mosquito Cells and Extends Viral Replication *Viruses* **9(9)**:262.

Theodosakis *et al.* (2017) p90RSK Blockade Inhibits Dual BRAF and MEK Inhibitor-Resistant Melanoma by Targeting Protein Synthesis. *J. Invest. Dermatol.* **137(10)**:2187.

Arimoto-Matsuzaki *et al.* (2016) TIA1 oxidation inhibits stress granule assembly and sensitizes cells to stress-induced apoptosis. *Nat. Commun.* **7**:10252.

Cinti *et al.* (2016) HIV-1 Gag Blocks Selenite-Induced Stress Granule Assembly by Altering the mRNA Cap-Binding Complex *mBio* **7(2)**:e00329-16.

Timpano *et al.* (2016) Analysis of Cap-binding Proteins in Human Cells Exposed to Physiological Oxygen Conditions. *J. Vis. Exp.* doi:10.3791.

Wahba *et al.* (2016) Polysome Profiling Links Translational Control to the Radioresponse of Glioblastoma Stem-like Cells. *Cancer Res.* **76(10)**:3078.

Hatakeyama *et al.* (2014) A Novel Functional Site in the PB2 Subunit of Influenza A Virus Essential for Acetyl-CoA Interaction, RNA Polymerase Activity, and Viral Replication. *J. of Biol. Chem.* **289**:24980.

Igreja *et al.* (2014) 4E-BPs require non-canonical 4E-binding motifs and a lateral surface of eIF4E to repress translation *Nat. Commun.* **5**:4790.

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

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