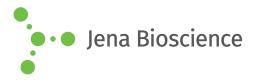
DATA SHEET

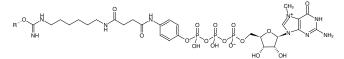




Immobilized γ-Aminophenyl-m⁷GTP (C₁₀-spacer)

y-Aminophenyl-m⁷GTP (C₁₀-spacer)-Agarose

Amount
1 ml
5 ml



Structural formula of Immobilized γ-Aminophenyl-m⁷GTP (C₁₀-spacer)

	Agarose characteristics
Bead/Particle size	45-165 μm
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 µmol - 7 µmol m⁷GTP/ml gel

Storage buffer: 20% Ethanol

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.

