



## Immobilized 2'/3'-EDA-m<sup>7</sup>GTP

2'/3'-EDA-7-Methyl-guanosine 5'-triphosphate (2'/3'-EDA-m<sup>7</sup>GTP) immobilized on Agarose  
2'/3'-EDA-m<sup>7</sup>GTP-Agarose

| Cat. No. | Amount |
|----------|--------|
| AC-142S  | 1 ml   |
| AC-142L  | 5 ml   |

**For research use only!**

**Shipping:** shipped at 4 °C

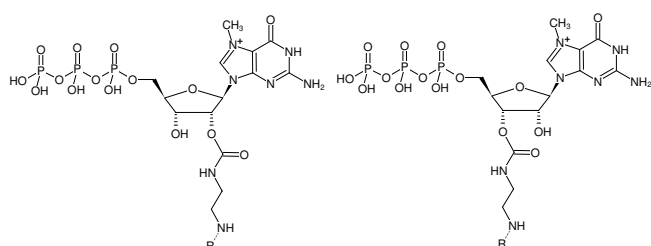
**Storage Conditions:** store at 4 °C

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

**Shelf Life:** 12 months after date of delivery

**Degree of substitution:** 5 µmol - 7 µmol 2'/3'-EDA-m<sup>7</sup>GTP/ml gel

**Storage buffer:** 20% Ethanol



Structural formula of Immobilized 2'/3'-EDA-m<sup>7</sup>GTP

|                              | 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<br><b>Not stable in organic solvents!</b> |
| Sterilization                | Not autoclavable!  |

R= Agarose

### Selected References:

Zhu *et al.* (2019) Targeting KPNB1 overcomes TRAIL resistance by regulating DR5, Mcl-1 and FLIP in glioblastoma cells. *Cell Death Dis.* **10** (2):118.

Zhu *et al.* (2019) XPO1 inhibitor KPT-330 synergizes with Bcl-xL inhibitor to induce cancer cell apoptosis by perturbing rRNA processing and Mcl-1 protein synthesis. *Cell Death Dis.* **10** (6):395.

William *et al.* (2018) eIF4E-Binding Proteins 1 and 2 Limit Macrophage Anti-Inflammatory Responses through Translational Repression of IL-10 and Cyclooxygenase-2. *J. Immunol.* **doi:10.4049.**

Posternak *et al.* (2017) MYC Mediates mRNA Cap Methylation of Canonical Wnt/β-catenin Signaling Transcripts by Recruiting CDK7 and RNA Methyltransferase. *Mol. Cancer Res.* **15** (2):213.

Xu *et al.* (2016) Improved transcription and translation with L-leucine stimulation of mTORC1 in Roberts syndrome. *BMC Genomics* **17**:25.

Liberman *et al.* (2015) DAP5 associates with eIF2β and eIF4AI to promote Internal Ribosome Entry Site driven translation. *Nucleic Acids Res.* **43** (7):3764.