



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

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

Cat. No.	Amount
AC-143S	1 ml
AC-143L	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. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

**Shelf Life:** 12 months

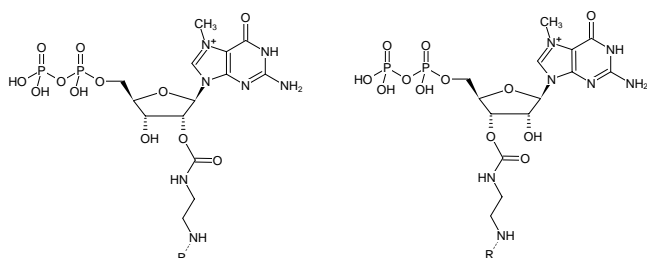
**Degree of substitution:** 5 μmol 2'/3'-EDA-m<sup>7</sup>-GDP/ml gel

**Storage buffer:** 20% Ethanol

**Please note:** For the purification of eukaryotic mRNA cap-binding proteins we recommend AC-155.

### Selected References:

Liu *et al.* (2015) RACK1-mediated translation control promotes liver fibrogenesis. *Biochem. Biophys. Res. Commun.* **463** (3):255.



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

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