



Immobilized 2'/3'-EDA-m⁷GDP

2'/3'-EDA-7-Methyl-guanosine 5'-diphosphate (2'/3'-EDA-m⁷GDP) immobilized on Agarose
2'/3'-EDA-m⁷GDP-Agarose

Cat. No.	Amount
AC-143L	5 ml
AC-143S	1 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

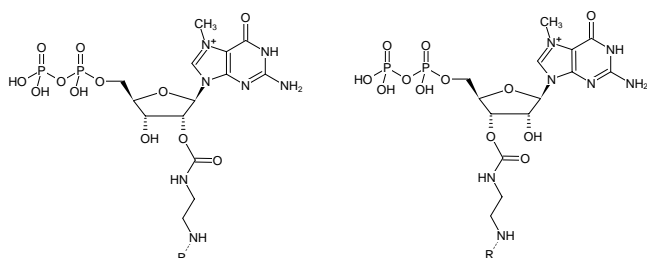
Degree of substitution: 5 μmol 2'/3'-EDA-m⁷-GDP/ml gel

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-m⁷GDP

	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!