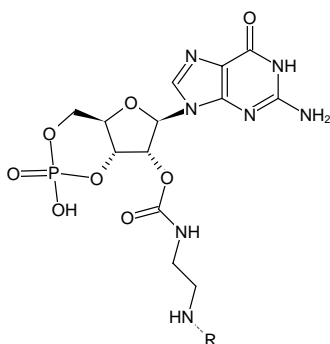




Immobilized 2'-EDA-cGMP

2'-O-(2-Aminoethyl-carbamoyl)-guanosine-3',5'-cyclic-monophosphate (EDA-cGMP) immobilized on Agarose.
2'-EDA-cGMP-Agarose

Cat. No.	Amount
AC-148S	1 ml
AC-148L	5 ml



Structural formula of Immobilized 2'-EDA-cGMP

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 EDA-cGMP/ml gel

Storage buffer: 20% Ethanol

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