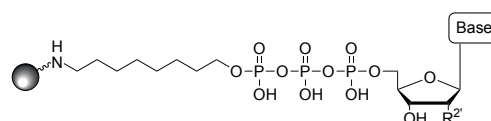


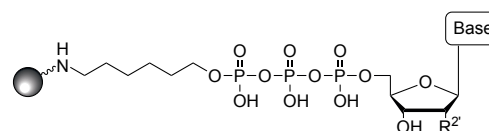
Immobilized Nucleotides: The perfect match!

γ -Phosphate-Amino-octyl-Linkage

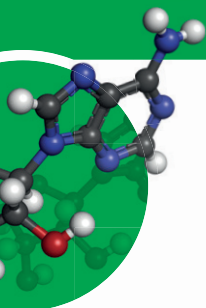


Affinity resin	Cat-No.	Amount
Immobilized γ -Amino-octyl-ATP	AC-105S	1 ml
	AC-105L	5 ml
Immobilized γ -Amino-octyl-GTP ^[1]	AC-106S	1 ml
	AC-106L	5 ml
Immobilized γ -Amino-octyl-dATP	AC-111S	1 ml
	AC-111L	5 ml
Immobilized γ -Amino-octyl-dGTP	AC-112S	1 ml
	AC-112L	5 ml

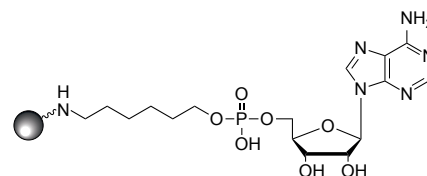
γ -Phosphate-Amino-hexyl-Linkage



Affinity resin	Cat-No.	Amount
Immobilized γ -Amino-hexyl-ATP	AC-116S	1 ml
	AC-116L	5 ml
Immobilized γ -Amino-hexyl-GTP ^[2]	AC-117S	1 ml
	AC-117L	5 ml
Immobilized γ -Amino-hexyl-dATP ^[3]	AC-122S	1 ml
	AC-122L	5 ml
Immobilized γ -Amino-hexyl-dGTP	AC-123S	1 ml
	AC-123L	5 ml
Immobilized γ -Amino-hexyl-m ⁷ GTP ^[4]	AC-141S	1 ml
	AC-141L	5 ml

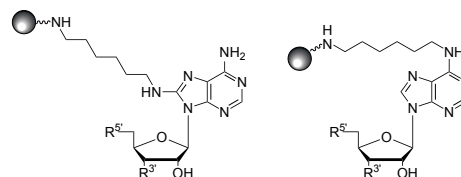


α -Phosphate-Amino-hexyl-Linkage



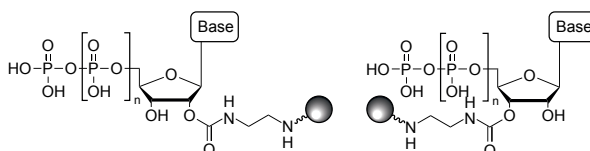
Affinity resin	Cat-No.	Amount
Immobilized α -Amino-hexyl-AMP	AC-158S	1 ml
	AC-158L	5 ml

Base Linkage

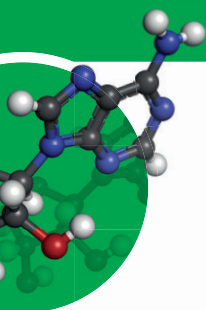


Affinity resin	Cat-No.	Amount
Immobilized 8-Amino-hexyl-ATP	AC-127S	1 ml
	AC-127L	5 ml
Immobilized 8-Amino-hexyl-cAMP	AC-146S	1 ml
	AC-146L	5 ml
Immobilized 8-Amino-hexyl-AMP	AC-156S	1 ml
	AC-156L	5 ml
Immobilized N ⁶ -(6-Amino)hexyl-ATP ^[5]	AC-129S	1 ml
	AC-129S	5 ml
Immobilized N ⁶ -(6-Amino)hexyl-AMP (identical to 5' AMP Sepharose 4B) ^[6]	AC-145S	1 ml
	AC-145L	5 ml

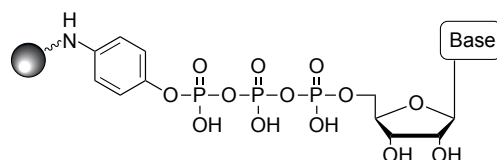
2'/3'-EDA-Linkage



Affinity resin	Cat-No.	Amount
Immobilized 2'/3'-EDA-ATP ^[5]	AC-131S	1 ml
	AC-131L	5 ml
Immobilized 2'/3'-EDA-GTP	AC-132S	1 ml
	AC-132L	5 ml
Immobilized 2'/3'-EDA-m ⁷ GTP ^[7,8]	AC-142S	1 ml
	AC-142L	5 ml
Immobilized 2'/3'-EDA-m ⁷ GDP	AC-143S	1 ml
	AC-143L	5 ml
Immobilized 2'/3'-EDA-AMP	AC-157S	1 ml
	AC-157L	5 ml
Immobilized 2'-EDA-cAMP	AC-147S	1 ml
	AC-147L	5 ml
Immobilized 2'-EDA-cGMP	AC-148S	1 ml
	AC-148L	5 ml



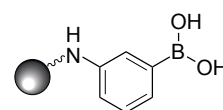
γ-Aminophenyl-Linkage



Affinity resin	Cat-No.	Amount
Immobilized γ-Aminophenyl-ATP (C ₁₀ -spacer) ^[9,10]	AC-101S	1 ml
	AC-101L	5 ml
Immobilized γ-Aminophenyl-ATP (no spacer)	AC-102S	1 ml
	AC-102L	5 ml
Immobilized γ-Aminophenyl-m ⁷ GTP (C ₁₀ -spacer) ^[11-13]	AC-155S	1 ml
	AC-155L	5 ml

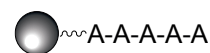
General purpose resin

for the enrichment and isolation of 1,2-cis-diols such as ribonucleotides, cofactors and carbohydrates



Affinity resin	Cat-No.	Amount
Immobilized m-Aminophenylboronic acid	AC-160	5 ml

Isolation of Poly(A)-binding Proteins

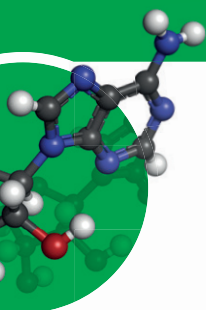


Affinity resin	Cat-No.	Amount
Poly(A)-Agarose	AK-159S	1 ml
	AK-159L	5 ml

Affinity Chromatography Kits for Adenosine Nucleotide binding Proteins

fast and easy purification

Affinity resin	Cat-No.	Amount
AMP Affinity Test Kit	AK-107	1 Kit
AMP Affinity Test Kit ^[14,15]	AK-102	1 Kit
AMP AffiPur Test Kit I ^[16,17]	AK-103S	1 Kit
	AK-103L	1 Kit
AMP AffiPur Test Kit II	AK-104S	1 Kit
	AK-104L	1 Kit
AMP AffiPur Test Kit III	AK-105S	1 Kit
	AK-105L	1 Kit
AMP AffiPur Test Kit IV	AK-106S	1 Kit
	AK-106L	1 Kit



Affinity Kit Components

all kit components are available separately

Affinity resin	Cat-No.	Amount
ATP Binding Buffer (5×)	AK-102B-S	15 ml
	AK-102B-L	75 ml
ATP WASH Buffer (5×)	AK-102W-S	10 ml
	AK-102W-L	50 ml
ATP Elution Buffer (5×)	AK-102E-S	2 ml
	AK-102E-L	10 ml
Protease Inhibitor Mix (100×)	AK-102I-S	0,5 ml
	AK-102I-L	5 × 5 ml
Sodium Orthovanadate (100×)	AK-102V-S	100 µl
	AK-102V-L	500 µl
PBS Tablets	AK-102P-S	10 tablets
	AK-102P-L	15 tablets

Base Linkage

- [1] Yun *et al.* (2015) An early endosome regulator, Rab5b, is an LRRK2 kinase substrate. *J. Biochem.* **157** (6):485.
- [2] Deng *et al.* (2019) Ubiquitination of Rheb governs growth factor-induced mTORC1 activation. *Cell Res.* **29** (2):136.
- [3] Domkin *et al.* (2014) Phosphines are ribonucleotide reductase reductants that act via C-terminal cysteines similar to thioredoxins and glutaredoxins. *Sci. Rep.* **4**:5539.
- [4] Landon *et al.* (2014) MNKS act as a regulatory switch for eIF4E1 and eIF4E3 driven mRNA translation in DLBCL. *Nat. Commun.* **5**:5413.
- [5] Park *et al.* (2015) PUGNAc induces protein ubiquitination in C2C12 myotube cells. *Cell Biochem. Funct.* **33** (8):525.
- [6] Yegutkin *et al.* (2014) Consequences of the lack of CD73 and prostatic acid phosphatase in the lymphoid organs. *Mediators Inflamm.* **2014**:485743.
- [7] Liberman *et al.* (2015) DAP5 associates with eIF2β and eIF4AI to promote Internal Ribosome Entry Site driven translation. *Nucleic Acids Res.* **43** (7):3764.
- [8] 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.* **200** (12):4102.
- [9] Nariai *et al.* (2015) Disruption of Heat Shock Protein 90 (Hsp90)-Protein Kinase Cδ (PKCδ) Interaction by (-)-Maackiain Suppresses Histamine H1 Receptor Gene Transcription in HeLa Cells. *J. Biol. Chem.* **290** (45):27393.
- [10] Teixeira-Duarte *et al.* (2019) Activation of a nucleotide-dependent RCK domain requires binding of a cation cofactor to a conserved site. *eLife* **8**:e50661.
- [11] Choe *et al.* (2018) mRNA circularization by METTL3-eIF3h enhances translation and promotes oncogenesis. *Nature* **561** (7724):556.
- [12] Grüner *et al.* (2016) The Structures of eIF4E-eIF4G Complexes Reveal an Extended Interface to Regulate Translation Initiation. *Mol. Cell.* **64** (3):467.
- [13] Goldsmith *et al.* (2020) Ribosome profiling reveals a functional role for autophagy in mRNA translational control. *Commun. Biol.* **3**:388.
- [14] Jain *et al.* (2020) Development of novel anti-malarial from structurally diverse library of molecules, targeting plant-like Calcium Dependent Protein Kinase 1, a multistage growth regulator of *P. falciparum*. *Biochem. J.* **477** (10):1951.
- [15] Nganon *et al.* (2019) *Plasmodium* pseudo-Tyrosine Kinase-like binds PP1 and SERA5 and is exported to host erythrocytes. *Sci. Rep.* **9**:8120.
- [16] Ding *et al.* (2015) The Q Motif Is Involved in DNA Binding but Not ATP Binding in ChIR1 Helicase. *PLoS One* **10** (10):e0140755.
- [17] Li *et al.* (2014) Synphilin-1 Binds ATP and Regulates Intracellular Energy Status. *PLoS One* **9** (12):e115233.

