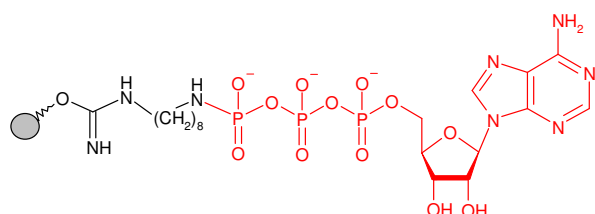


## $\gamma$ -Amino-octyl-ATP - Sepharose<sup>®</sup>

Cat. No.	Amount
AC-105S	1 ml (bulk material)
AC-105L	5 ml (bulk material)
AC-105C	1 ml (syringe column)
AC-105SC	0.2 ml (scr. column)
AC-105MP	1 ml (MPLC column)



Adenosine triphosphate (ATP) immobilized on Agarose, suitable for purification of ATP-binding proteins.

**Degree of substitution:** 5  $\mu$ mol ATP/ml Gel

**Linker:** 10 atoms

**Attachment:**  $\gamma$ -phosphate

**Storage buffer:** 100 mM Tris, pH 7.5, 50 % Glycerol, 0.02 % Thimerosal

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

Long term storage at 4°C.

**Expiry:** 12 month

<sup>®</sup> Sepharose is a trademark of Amersham-Biosciences.

### Selected References:

Suto *et. al.* (1998) Synthesis of  $\gamma$ -phosphate linked nucleoside affinity chromatography resins for protein purification, including ribonucleoside triphosphate reductase. *Nucleosides and nucleotides*, **17(8)**:1453