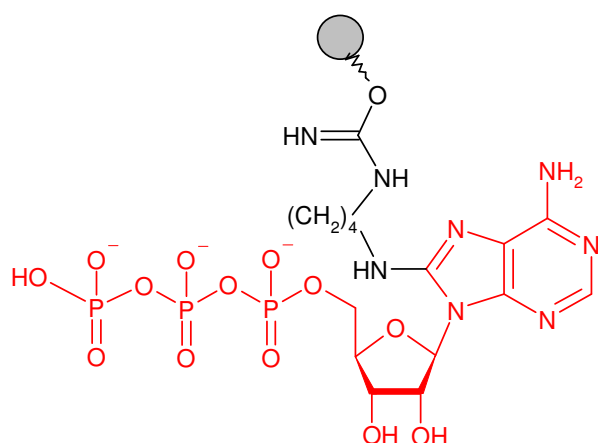


## 8-Amino-butyl-ATP - Sepharose®

Cat. No.	Amount
AC-128S	1 ml (bulk material)
AC-128L	5 ml (bulk material)
AC-128C	1 ml (syringe column)
AC-128SC	0.2 ml (scr. column)
AC-128MP	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:** 6 atoms

**Attachment:** C8

**Storage buffer:** 50% glycerol (cont. 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

® Sepharose is a trademark of Amersham-Biosciences.