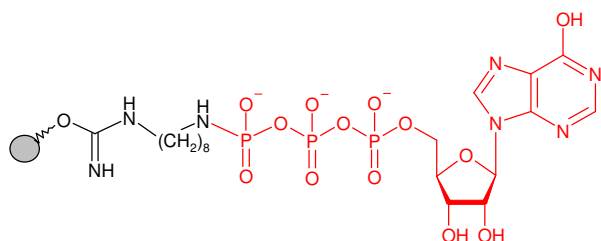


γ -Amino-octyl-ITP - Sepharose[®]

Cat. No.	Amount
AC-109S	1 ml (bulk material)
AC-109L	5 ml (bulk material)
AC-109C	1 ml (syringe column)
AC-109SC	0.2 ml (scr. column)
AC-109MP	1 ml (MPLC column)



Inosine 5'-triphosphate (ITP) immobilized on Agarose, suitable for purification of ITP-binding proteins.

Degree of substitution: 5 μ mol ITP/ml Gel

Linker: 10 atoms

Attachment: γ -phosphate

Storage buffer: 20% Ethanol

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.

Selected References:

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