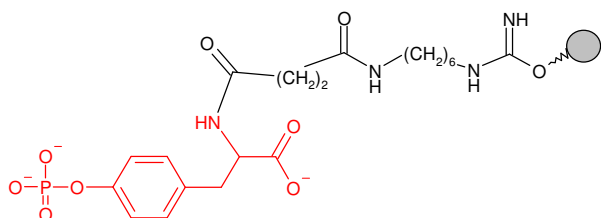


O-Phospho-L-Tyrosine - Sepharose® (C₁₀-spacer)

Cat. No.	Amount
AC-103S	1 ml (bulk material)
AC-103L	5 ml (bulk material)
AC-103C	1 ml (syringe column)
AC-103SC	0.2 ml (scr. column)
AC-103MP	1 ml (MPLC column)



O-Phospho-L-Tyrosine immobilized on Agarose, suitable for purification of proteins containing a SH2-domain.

Degree of substitution: 25 μmol O-Phospho-L-Tyrosine/ ml Gel

Linker: C₁₀

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:

Fry *et al.* (1992) Purification and characterization of a phosphatidyl 3-kinase complex from bovine brain by using phosphopeptide affinity columns. *Biochem. J.* **288**(2): 383.