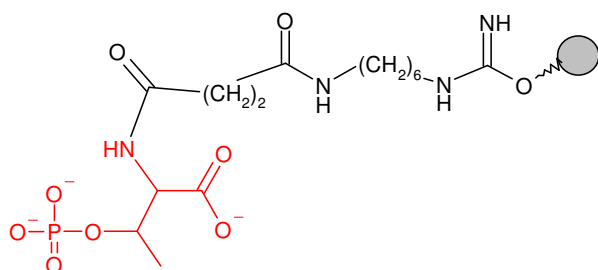


O-Phospho-L-Threonine - Sepharose® (C₁₀-spacer)

Cat. No.	Amount
AC-139S	1 ml (bulk material)
AC-139L	5 ml (bulk material)
AC-139C	1 ml (syringe column)
AC-139SC	0.2 ml (scr. column)
AC-139MP	1 ml (MPLC column)



O-Phospho-L-Threonine immobilized on Agarose, suitable for purification of protein threonine phosphatases.

Degree of substitution: 25 μmol O-Phospho-L-Threonine/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.