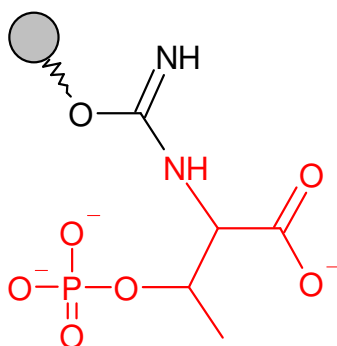


O-Phospho-L-Threonine - Sepharose®

(no spacer)

Cat. No.	Amount
AC-140S	1 ml (bulk material)
AC-140L	5 ml (bulk material)
AC-140C	1 ml (syringe column)
AC-140SC	0.2 ml (scr. column)
AC-140MP	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: none

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.