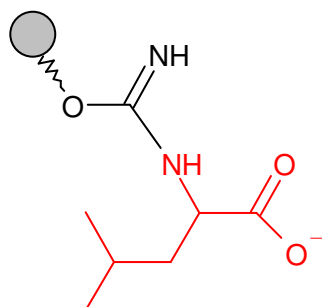


L-Leucine - Sepharose®

(no spacer)

Cat. No.	Amount
AC-151S	1 ml (bulk material)
AC-151L	5 ml (bulk material)
AC-151C	1 ml (syringe column)
AC-151SC	0.2 ml (scr. column)
AC-151MP	1 ml (MPLC column)



(S)-2-Amino-4-methylpentanoic acid, L-Leucine immobilized on Agarose

Degree of substitution: 5 μmol L-Leucine/ml Gel 5 μmol 2'/3'-EDA-m7-GDP/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.