

Separation of di-peptides

HPTLC Cellulose ohne F

Plate	HPTLC pre-coated plate cellulose, without indicator (P/N 1. 05651, alt. P/N 1. 16092)
Eluent	Ethanol/water 80/20 (v/v) with addition of 0,05 mol/l lithium chloride
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Alg-Asp 2. Ala-Gly 3. Ala-Ala
Detection	Spray reagent: Ninhydrin Heating up to 110°C for 5min.

