

Separation of di-peptides

HPTLC Silica gel 60 RP-2 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s (1. 13726)
Eluent	Acetonitrile + buffer solution pH 3.00 50ml+5ml with addition of 0,05 mol/l lithium chloride
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Ala-Gly 2. Ala-Ala 3. Asp-Leu
Detection	Spray reagent: Ninhydrin Heating up to 110°C for 10 min.

