

Separation of penta-peptides

TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N 1.15389, alt. P/N 1.15423)
Eluent	Acetonitrile/ethanol/IN hydrochloric acid, 85/5/10 (v/v/v) with addition of 0,1 mol/l tetramethylammonium chloride
Migration distance	12 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Glu-Thr-Tyr-Ser-Lys 2. Ala-Ala-Ala-Ala-Ala 3. Gly-Gly-Asn-Asp-Ala
Application volume	700 nl
Detection	Spray reagent: Ninhydrin Heating up to 120°C for 10 min.

