

Separation of penta-peptides

TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N: 1.15389, Alternative P/N: 1.15423)
Eluent	Acetonitrile/1N hydrochloric acid 90/10 (v/v) with addition of 0,1 mol/l tetramethylammonium chloride
Migration Distance	8 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 4. Ala-Ala-Tyr-Ala-Ala 5. Leu-Trp-Met-Arg-Phe (all 0,2 %)
Application volume	700 nl
Detection	Spray reagent: ninhydrin heating up to 120°C for 10 min.

