

Separation of tetra-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/water 40/60 (v/v) with addition of 0,1 mol/l lithium chloride
Migration distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Phe-Phe-Phe-Phe 2. Val-Val-Val-Val 3. Gly-Gly-Glu-Ala 4. Ala-Ala-Ala-Ala (all 0,4%)
Application volume	800 nl
Detection	Spray reagent: Ninhydrin Heating up to 120°C for 10 min.

