

Separation of tri-peptides

HPTLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-18 F 254s (1. 13725)
Eluent	Acetonitrile/buffer solution pH 3.00 90/10 (v/v) with addition of 0,1 mol/l lithium chloride
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Ala-Ala-Ala 2. Leu-Gly-Gly 3. Leu-Leu-Leu
Detection	Spray reagent: Ninhydrin Heating up to 110°C for 20 min.

