

# Separation of Fmoc-peptides

## HPTLC Silica gel 60 RP-18 WF254s

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Acetonitrile/water 50/50 (v/v) with addition of 0,1 mol/l lithium chloride
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Fmoc-Leu-Leu-Leu 2. Fmoc-Leu-Leu 3. Fmoc-Leucine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

