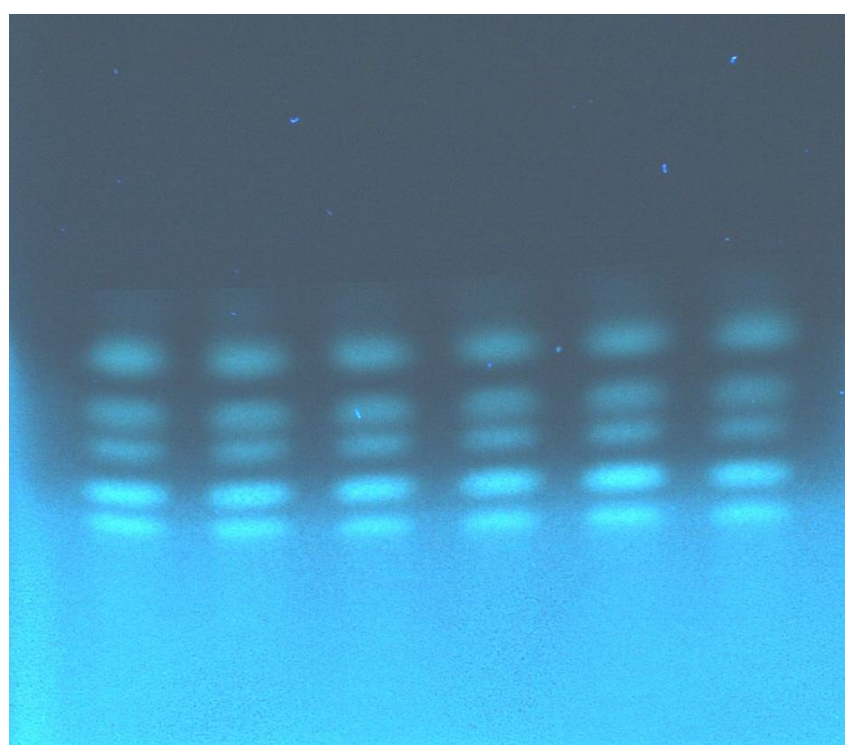


Separation of peptides

Plate	HPTLC Silica gel 60 RP-8 F254s (P/N 1. 13725)
Mobile phase (v/v)	methanol/water (65/35) + 0,1mol/l ammonium
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Detection	0,02% Fluorescamin solution in acetone spray 20s and dry 10min at roomtemperature Detection UV @ 366nm



← Gly-Tyr
 ← Val-Tyr-Val
 ← Methionine enkephalin
 ← Leucine enkephalin
 ← H-YGSG-pT-AK-NH3

Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent
1 to 6 peptide mix	H-YGSG-pT-AK-NH2	0,9	water/ACN 85/15
	leucine enkephalin	1,0	
	methionine enkephalin	1,0	
	Val-Try-Val	1,5	
	Gly-Try	1,2	