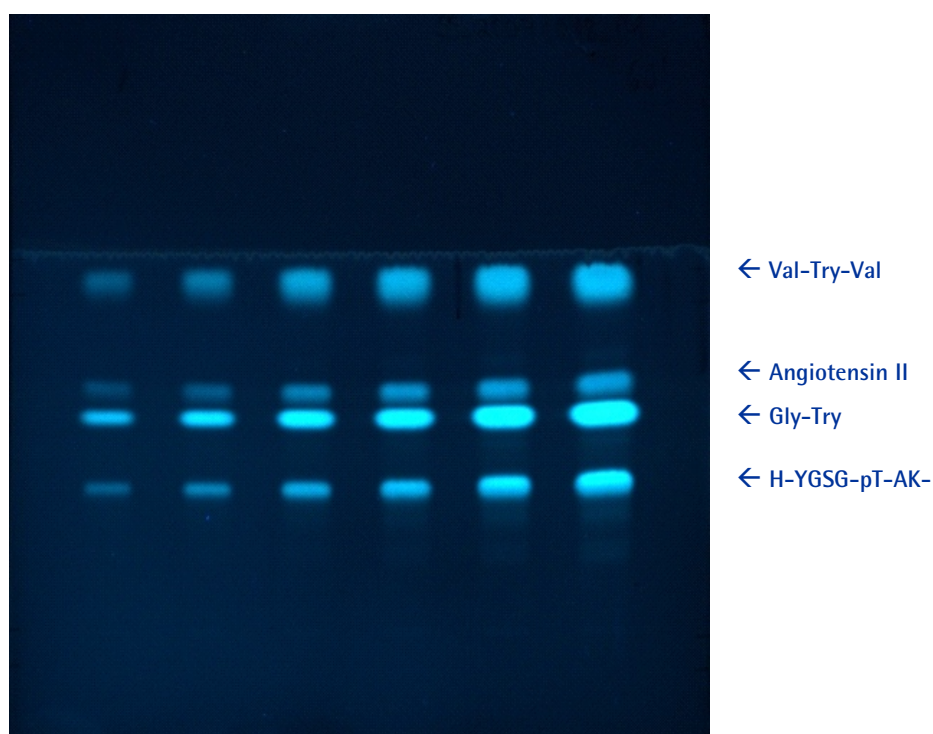


Separation of peptides on HPTLC Cellulose

Plate	ProteoChrom® HPTLC Cellulose (P/N 1. 05651)
Mobile phase (v/v)	2-butanol/pyridin/acetic acid/water (15/10/3/12)
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



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	Application volume µl
1 to 6	peptide mix			
	H-YGSG-pT-AK-NH ₂	1,0	water/ACN 85/15	0,1
	Gly-Try	1,0		0,5
	Val-Try-Val	1,0		1,0
	angiotensin II	1,2		3,0
				5,0
				7,0