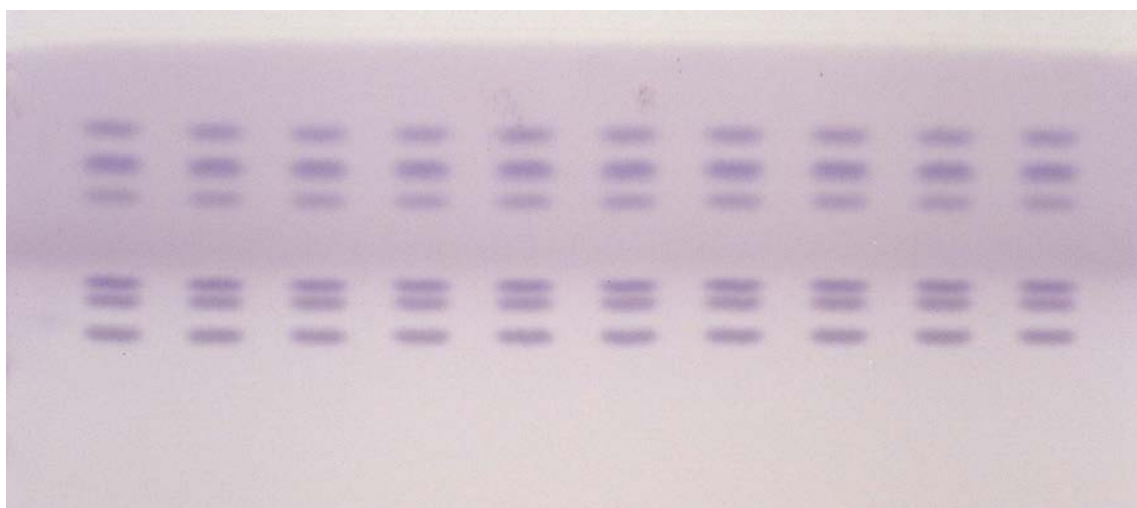


Separation of amino acids on Cellulose

Plate	TLC Cellulose 20x10cm (P/N 1. 05716, alt P/N 1. 05730)
Mobile phase (v/v)	1. 1-butanol/acetone/glacial acetic acid/water 2. 1-butanol/acetone/glacial acetic acid/water + 1ml Ninhydrin solution Ninhydrin solution: 7g Ninhydrin in 100ml acetone/1-
Migration distance	Each 7 cm
Application volume	2 µL
Chamber	Normal chamber without chamber saturation
Detection	Heat 1-2min at 110°C, Detection white light



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf	
1 to 10	Amino- acid- mix	L-serine	0,15	water/1-propanol 4/6	26
		L-threonine	0,15	water/1-propanol 4/6	34
		L-alanine	0,1	water/1-propanol 4/6	38
		L-methionine	0,15	water/1-propanol 4/6	58
		L-phenylalanine	0,25	water/1-propanol 4/6	67
		L-leucine	0,10	water/1-propanol 4/6	75