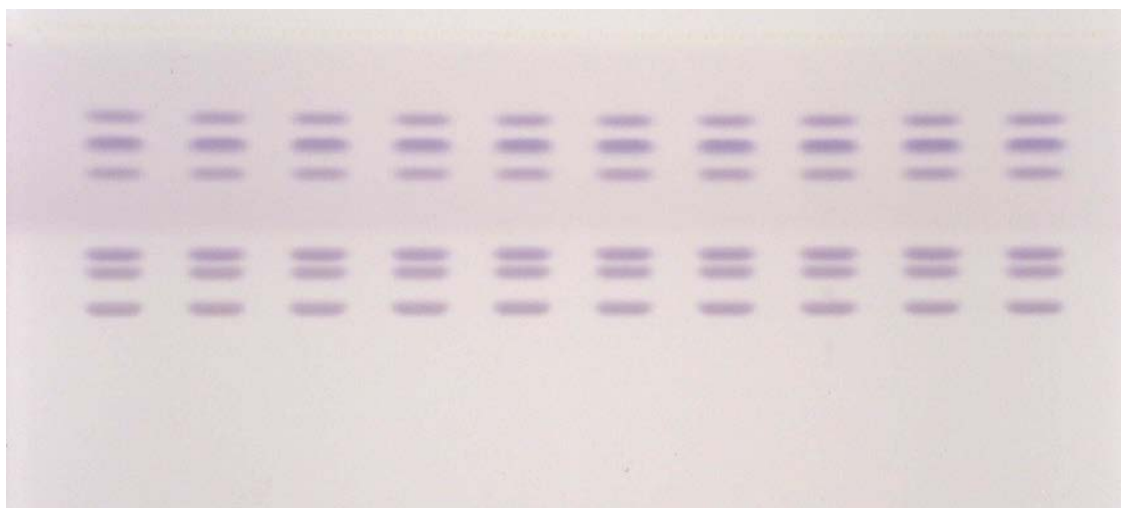


Separation of amino acids on Cellulose

Plate	HPTLC Cellulose 20x10cm (P/N 1. 05786, alt P/N 1. 05787)
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-Serin	0,15	water/1-propanol 4/6	31
		L-Threonin	0,15	water/1-propanol 4/6	40
		L-Alanin	0,1	water/1-propanol 4/6	45
		L-Methionin	0,15	water/1-propanol 4/6	65
		L-Phenylalanin	0,25	water/1-propanol 4/6	72
		L-Leucin	0,10	water/1-propanol 4/6	78