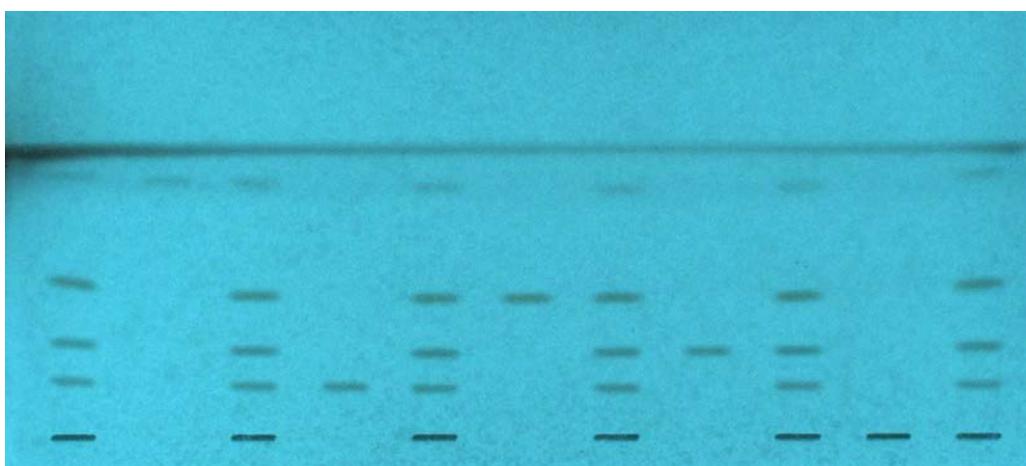


Separation of steroids Silica gel 60 CN

Plate	HPTLC Silica gel 60 CN F254s 20x10cm (P/N 1. 12571)
Mobile phase (v/v)	Petroleum benzine 40-60°C/acetone (80/20)
Migration distance	Each 5 cm
Application volume	1 µL
Chamber	Normal chamber without chamber saturation
Detection	20% perchloric acid in water spray 20s and heat 5min at 100°C, Detection UV @ 254nm



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	R _F
2	cholesterol	1,0	methanol	0
4	hydrocortisone	1,0	methanol	16
6	methyl testosterone	1,0	methanol	29
8	reichstein's S	1,0	methanol	48
10	trimipramine	1,0	methanol	86
1, 3, 5, 7, 9 and 11	steroid mixture	1,0	methanol	-