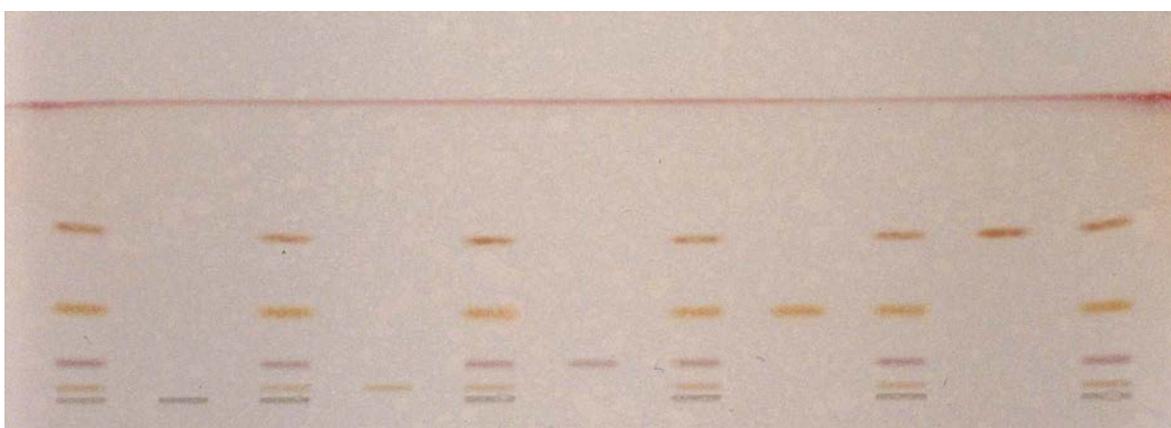


## Separation of steroids Silica gel 60 DIOL

Plate	HPTLC Silica gel 60 DIOL F254s 20x10cm (P/N 1. 05636, alt P/N 1. 12668)
Mobile phase (v/v)	Petroleum benzine 40–60°C/acetone (80/20)
Migration distance	Each 5 cm
Application volume	1 $\mu$ L
Chamber	Normal chamber without chamber saturation
Detection	20% perchloric acid in water spray 20s and heat 5min at 100°C, Detection white light



### Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
2	trimipramine	1,0	methanol	0
4	hydrocortisone	1,0	methanol	4
6	reichstein's S	1,0	methanol	13
8	methyl testosterone	1,0	methanol	31
10	cholesterol	1,0	methanol	56
1, 3, 5, 7, 9 and 11	steroid mixture	1,0	methanol	-