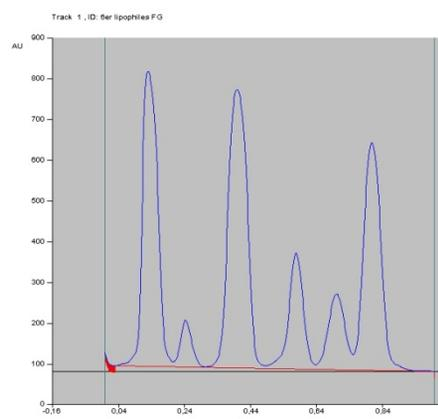
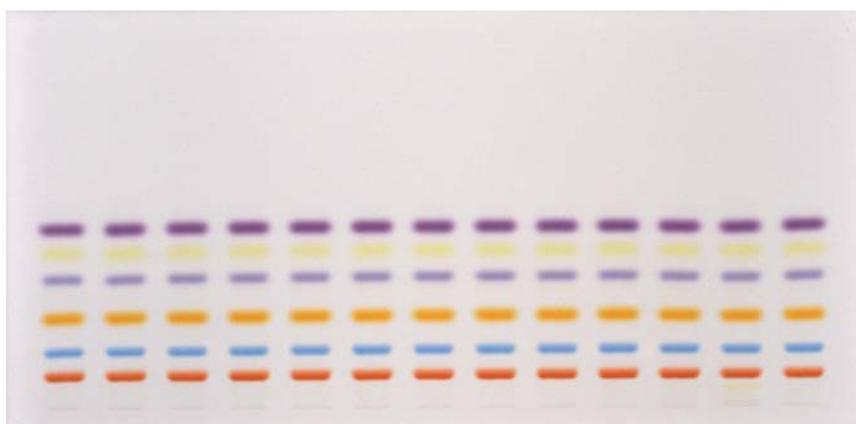


Separation of lipophilic dye mixture on Silica gel 60

Plate	HPTLC Silica gel 60 F254, 20x10cm (P/N 1. 05642)
Mobile phase (v/v)	Toluene
Migration distance	5 cm
Application volume	2 μ L
Chamber	Normal chamber without chamber saturation
Detection	Detection white light



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1 to 7	p-Methoxyazobenzol (52)	3,0	toluene	14
	Sudan II (79)	3,0	toluene	26
	Sudanviolett (89)	3,0	toluene	41
	Ceresviolett BRN (55)	1,0	toluene	59
				71
				82