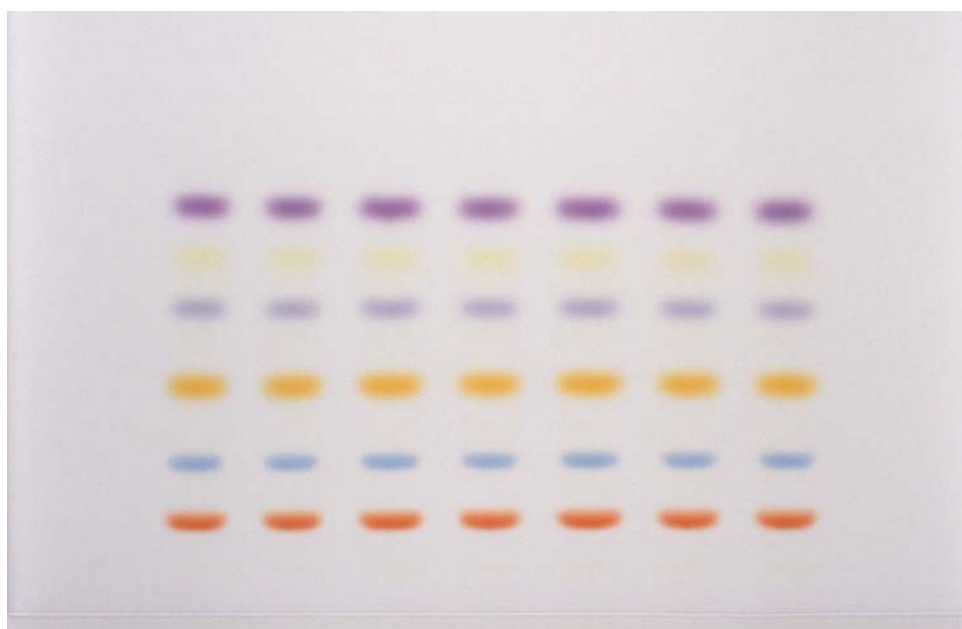


Separation of lipophilic dye mixture

Plate	TLC Silica gel 60 F254 with concentrating zone (P/N 1. 11798, alt P/N 1. 11846)
Mobile phase (v/v)	Toluene
Migration distance	10 cm
Application volume	2 µL
Chamber	Normal chamber without chamber saturation
Detection	Detection white light



Chromatographic data

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