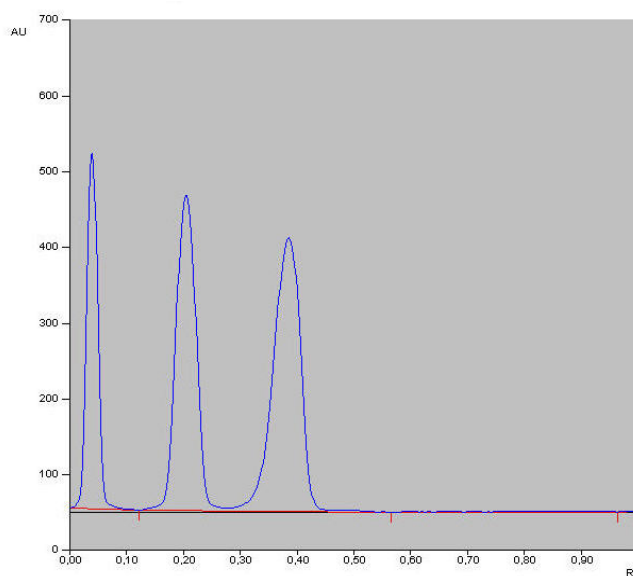


Separation of lipophilic dye mixture

Plate	HPTLC Silica gel 60 RP-18 F254s, 10x10cm (P/N 1. 13724, Alt P/N 1. 16225)
Mobile phase (v/v)	Acetonitrile/water(9/1)
Migration distance	5 cm
Application volume	500 nl
Chamber	Normal chamber without chamber saturation
Detection	Detection white light



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1 to 7	Fettrot 7B ()	1,0	acetone	4
	Ceresrot G ()	1,0	acetone	21
	Methylrot ()	0,2	acetone	40