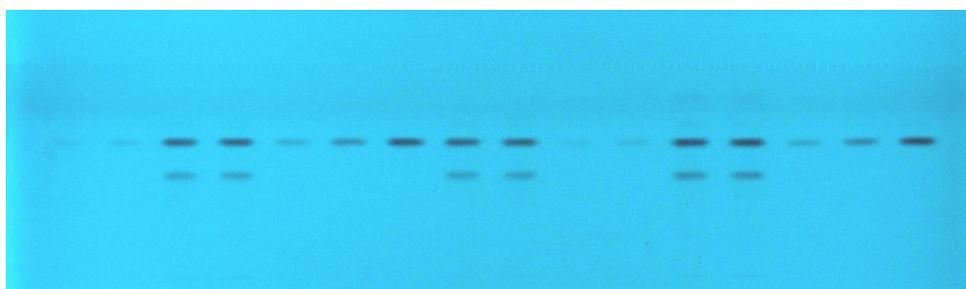
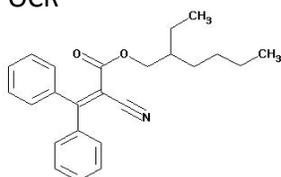


Separation of UV-Filter on Silica gel 60 RP-18

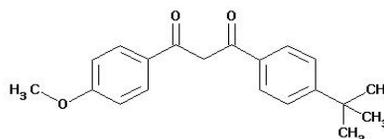
Plate	HPTLC Silica gel 60 RP-18 F254s 20x10cm (P/N 1. 16225, alt P/N 1. 13724)
Mobile phase (v/v)	Methanol/acetonitrile (90/10)
Migration distance	50 mm
Application volume	1 μ L
Chamber	Normal chamber without chamber saturation
Detection	Detection UV @ 254nm



Eusolex® OCR



Eusolex® 9020



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1 and 10	Eusolex® OCR	0,2	ethanol	64
2 and 11	Eusolex® OCR	0,4	ethanol	64
3 and 4	Eusolex® OCR and Eusolex® 9020, sample 1	-	ethanol	-
5 and 14	Eusolex® OCR	0,8	ethanol	64
6 and 15	Eusolex® OCR	1,6	ethanol	64
7 and 16	Eusolex® OCR	4,0	ethanol	64
8 and 9	Eusolex® OCR and Eusolex® 9020, sample 2	-	ethanol	-
12 and 13	Eusolex® OCR and Eusolex® 9020, sample 3	-	ethanol	-