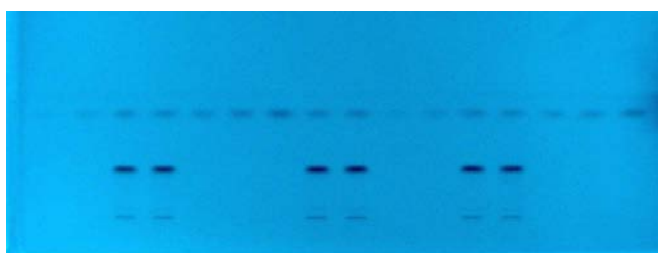
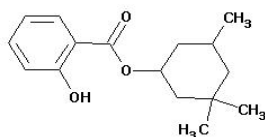


Separation of UV-Filter

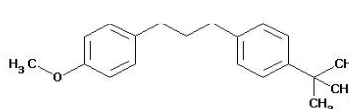
Plate	HPTLC Silica gel 60 F254s 20x10cm (P/N 1. 15696)
Mobile phase (v/v)	n-heptane/tetrahydrofuran (4/1)
Migration distance	50 mm
Application volume	1 µL
Chamber	Normal chamber without chamber saturation
Detection	Detection UV @ 254nm



Eusolex® HMS



Eusolex® 9020



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1 and 10	Eusolex® HMS	0,4	ethanol	84
2 and 11	Eusolex® HMS	1,0	ethanol	84
3 and 4	Eusolex® HMS, Eusolex 9020, Oxyxex ST and Ronacare AP sample 1	-	ethanol	-
5 and 14	Eusolex® HMS	1,5	ethanol	84
6 and 15	Eusolex® HMS	2,0	ethanol	84
7 and 16	Eusolex® HMS	3,0	ethanol	84
8 and 9	Eusolex® HMS, Eusolex 9020, Oxyxex ST and Ronacare AP sample 1	-	ethanol	-
12 and 13	Eusolex® HMS, Eusolex 9020, Oxyxex ST and Ronacare AP sample 1	-	ethanol	-