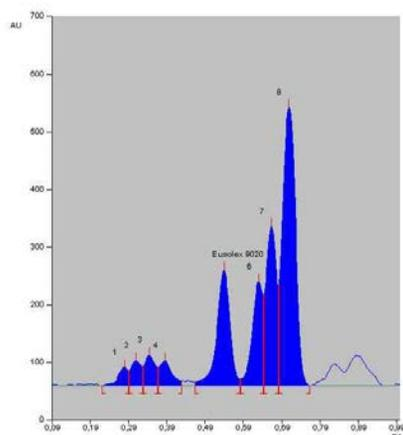
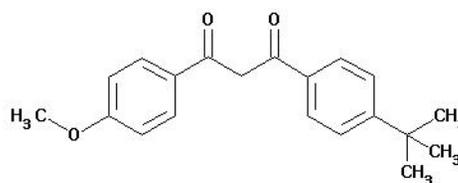


Separation of UV-Filter

Plate	HPTLC Silica gel 60 RP-18 F254s 20x10 cm (P/N 1. 16225, alt P/N 1. 13724)
Mobile phase (v/v)	Methanol
Application volume	1,5 µl
Migration distance	50 mm
Chamber	Normal chamber without chamber saturation
Detection	Detection UV @ 254nm



Eusolex® OCR



Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1, 6 and 11	Eusolex® 9020	0,2	ethanol	54
2, 7 and 12	Eusolex® 9020	0,5	ethanol	54
5, 10 and 15	Eusolex® 9020	0,8	ethanol	54
3, 4 and 8	Eusolex® OCR, Eusolex® 9020, Oxyhex ST, Ronacare AP, Finsolv TN	-	acetonitril	-
9, 13 and 14	Eusolex® OCR, Eusolex® 9020, Oxyhex ST, Ronacare AP, Finsolv TN	-	ethanol	-