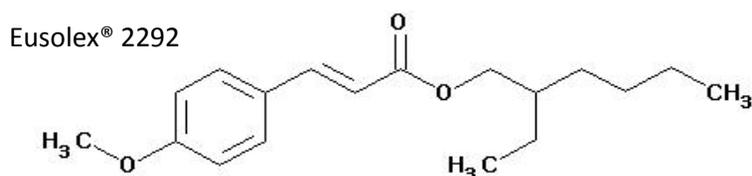
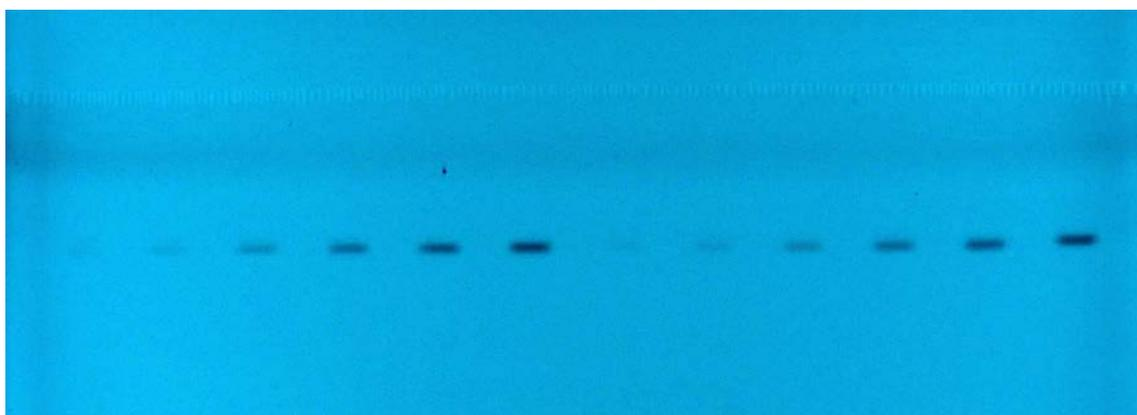


# Separation of UV-Filter

Plate	HPTLC Silica gel 60 RP-18 F254s 20x10 cm (P/N 1. 16225, alt P/N 1. 13724)
Mobile phase (v/v)	Methanol/ACN (90/10)
Migration distance	50 mm
Chamber	Normal chamber without chamber saturation
Detection	Detection UV @ 254nm



## Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf	Application volume $\mu$ l
1 and 10	Eusolex® 2292	2,0	methanol	46	0,1
2 and 11	Eusolex® 2292	2,0	methanol	46	0,2
3 and 4	Eusolex® 2292	2,0	methanol	46	0,5
5 and 14	Eusolex® 2292	2,0	methanol	46	1,0
6 and 15	Eusolex® 2292	2,0	methanol	46	1,5
7 and 16	Eusolex® 2292	2,0	methanol	46	2,0