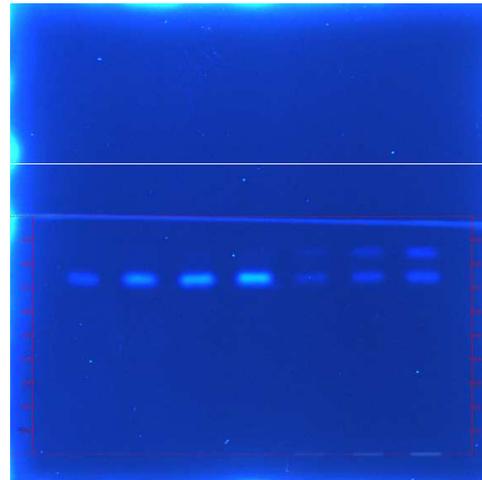
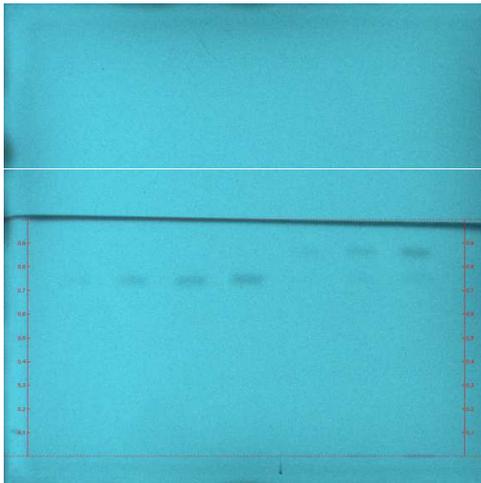




# Separation of a natural product from cosmetic matrix on HPTLC Silica gel 60 NH<sub>2</sub> F<sub>254</sub>S

## Chromatographic Conditions

Plate: HPTLC Silica gel 60 NH<sub>2</sub> F<sub>254</sub>S 10x10 cm  
 Article number: 1.15647.0001  
 Mobile Phase (v/v): Dichloremethan / Isopropanol 9:1  
 Migration distance: 5 cm  
 Chamber: Normal chamber without chamber saturation  
 Detection: UV 254 nm and UV 366 nm  
 Sample preparation: Standard was diluted in 10 ml isopropanol, sample was directly applied to the plate by CAMAG ATS4 6 mm bandwise.



## Chromatographic Data

Track No.	Compounds	Concentration (mg/ml)	Solvent	Application volume µl	hRF
1	Standard cos-active	1	Isopropanol	0.1	74
2	Standard cos-active	1	Isopropanol	0.2	74
3	Standard cos-active	1	Isopropanol	0.3	74
4	Standard cos-active	1	Isopropanol	0.4	74
5	Sample	~ 0.05		1	74 / 84
6	Sample	~ 0.05		2	74 / 84
7	Sample	~ 0.05		4	74 / 84