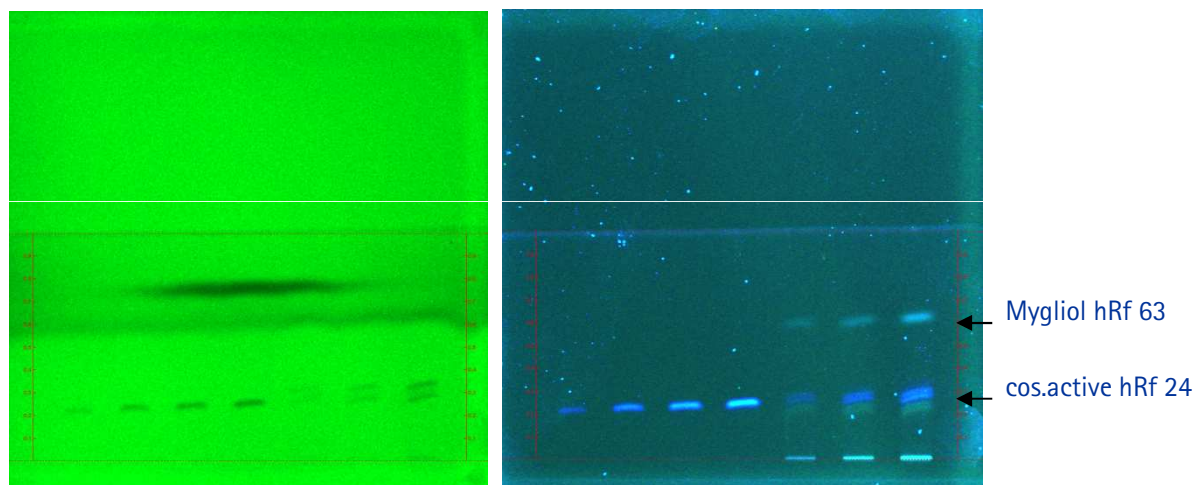




Determination of a cosmetic active from cosmetic oil on HPTLC Silica gel 60 F₂₅₄

Chromatographic Conditions

Plate: HPTLC Silica gel 60 F₂₅₄ 10x10 cm
 Article number: 1.05628.0001
 Mobile Phase (v/v): Dichloremethan
 Migration distance: 5 cm
 Chamber: Normal chamber without chamber saturation
 Detection: UV 254 nm, 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/10ml)	Solvent	Application volume µl	hRF
1	Cos-active- standard	1	Isopropanol	0.1	24
2	Cos-active- standard	1	Isopropanol	0.2	24
3	Cos-active- standard	1	Isopropanol	0.3	24
4	Cos-active- standard	1	Isopropanol	0.4	24
5	Cos-active sample	5 % in Mygliol		1	24/63
6	Cos-active sample	5 % in Mygliol		2	24/63
7	Cos-active sample	5 % in Mygliol		4	24/63