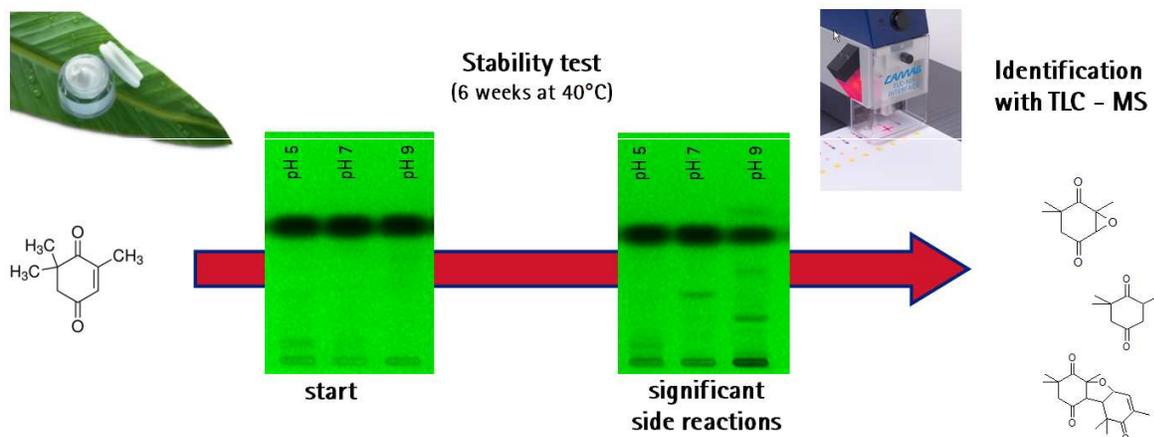


Determination of Oxoisophorone in cosmetic formulation on HPTLC Silica gel 60 F₂₅₄

Chromatographic Conditions

Plate:	HPTLC Silica gel 60 F ₂₅₄ 20x10 cm
Article number:	1.05649.0001
Mobile Phase (v/v):	Heptane / THF 8:2
Migration distance:	5 cm
Chamber:	Normal chamber without chamber saturation
Detection:	UV 254 nm, Identification of side products with TLC-MS Interface from CAMAG
Sample preparation:	1 g cosmetic formulation was stirred in a mixture of n-Heptane / THF 1:1 for 30 minutes and filtered over an 0.45 µm syringe filter. Application by CAMAG ATS4 6 mm bandwise.



Chromatographic Data

Track No.	Compounds	Concentration (mg/10ml)	Solvent	Application volume µl	hRF
1	Oxoisopherone in formulation pH 5	~10	Heptan/THF 1:1	2	56
2	Oxoisopherone in formulation pH 7	~10	Heptan/THF 1:1	2	56
3	Oxoisopherone in formulation pH 9	~10	Heptan/THF 1:1	2	56