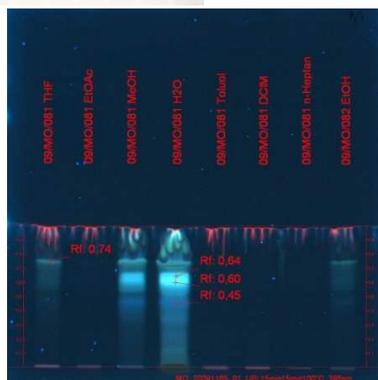
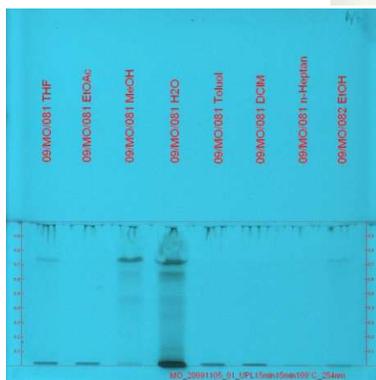


# Screening of plant extracts on HPTLC Silica gel 60 NH<sub>2</sub> F<sub>254</sub>S

## 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): Ethylacetat / Formic acid / Glacial acid / Water 100:11:11:27 + 0.1 % TFA  
 Migration distance: 5 cm  
 Chamber: Normal chamber without chamber saturation  
 Detection: UV 254 nm, UV 366 nm and white light  
 Sample preparation: 2 g plant material extracted for 1 h at 60°C under shaking with 50 ml solvent. Filtration of crude extract with an 0.45 µm syringe filter. Application by CAMAG ATS4 6 mm bandwise.



## Chromatographic Data

Track No.	Compounds	Concentration (mg/ml)	Solvent	Application volume µl
1	Extract	40	THF	2
2	Extract	40	Ethylacetate	2
3	Extract	40	Methanol	2
4	Extract	40	Water	2
5	Extract	40	Toluol	2
6	Extract	40	Methylechloride	2
7	Extract	40	n-Heptane	2
8	Extract	40	Methanol	2