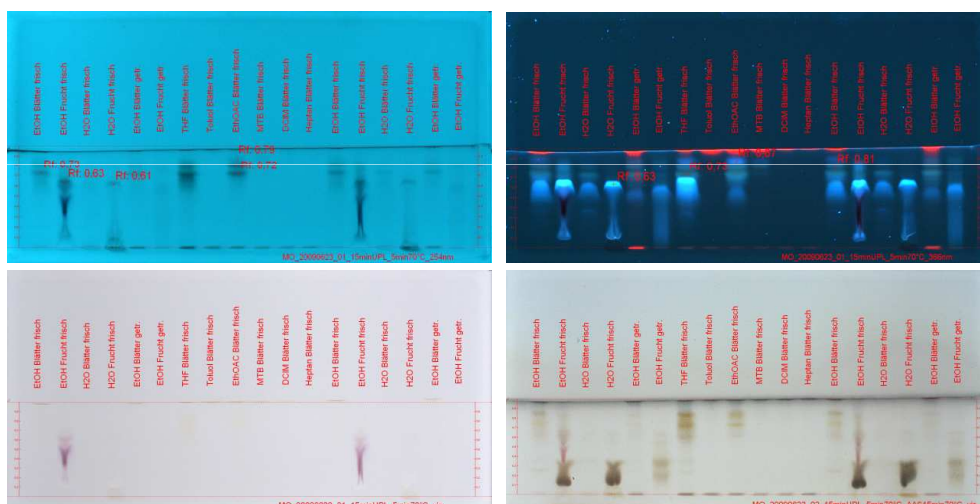




Detection of Compounds in amelanchir ovalis extract on HPTLC Silica gel 60 F₂₅₄S

Chromatographic Conditions

Plate: HPTLC Silica gel 60 F₂₅₄S 20x10 cm
 Article number: 1.15696.0001
 Mobile Phase (v/v): Ethylacetat / Formic acid / Glacial acid / Water 100:11:11:27 +0.1 % n-Heptane
 Migration distance: 5 cm
 Chamber: Normal chamber without chamber saturation
 Detection: UV 254 nm, UV 366 nm and white light, white light after derivatisation with Anisaldehyde- sulfuric acid-reagent
 Sample preparation: 1.9 g grounded leaves / 2,5 g smashed fruits were extracted for 30 h at 70°C with 10 ml solvent in an ultrasonic bath. Filtration of crude extract with an 0.45 µm syringe filter. Application by CAMAG ATS4 6 mm bandwise.



Chromatographic Data

Track No.	Compounds	Solvent	Application volume µl
1,2,5,6	Extract	Ethanol	2
3, 4	Extract	Water	2
7	Extract	Tetrahydrofuran	2
8	Extract	Toluol	2
9	Extract	Ethylacetate	2
10	Extract	MTB Ether	2
11	Extract	Dichlore methane	2
12	Extract	n-Heptane	2
13,14, 17,18	Extract	Ethanol	1
15, 16	Extract	Water	1