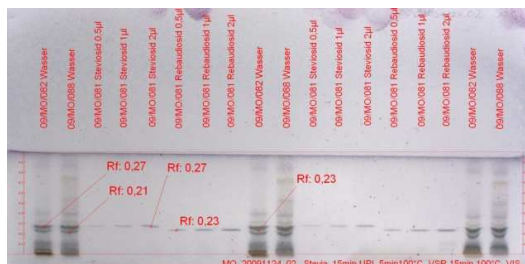
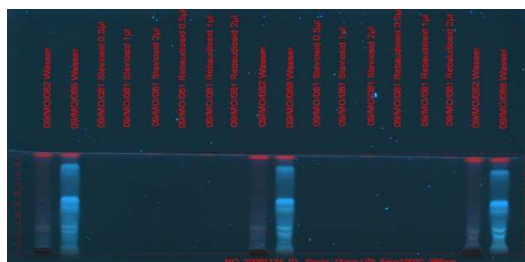




Detection of Stevioside and Rebaudioside in Stevia extract on HPTLC Silica gel 60 F₂₅₄S

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 % TFA
 Migration distance: 5 cm
 Chamber: Normal chamber without chamber saturation
 Detection: UV 254 nm, UV 366 nm, derivatised with Vanillin-sulfuric acid-reagent and Anisaldehyde-sulfuric acid-reagent detected with 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	hRF
1,9,17	Stevia extract 1	~0.04	Water	2	
2,10,18	Stevia extract 2	~0.04	Water	2	
3, 11	Stevioside standard	1	Water	0.5	27
4, 12	Stevioside standard	1	Water	1	27
5, 13	Stevioside standard	1	Water	2	27
6, 14	Rebaudioside standard	1	Water	0.5	23
7, 15	Rebaudioside standard	1	Water	1	23
8, 16	Rebaudioside standard	1	Water	2	23