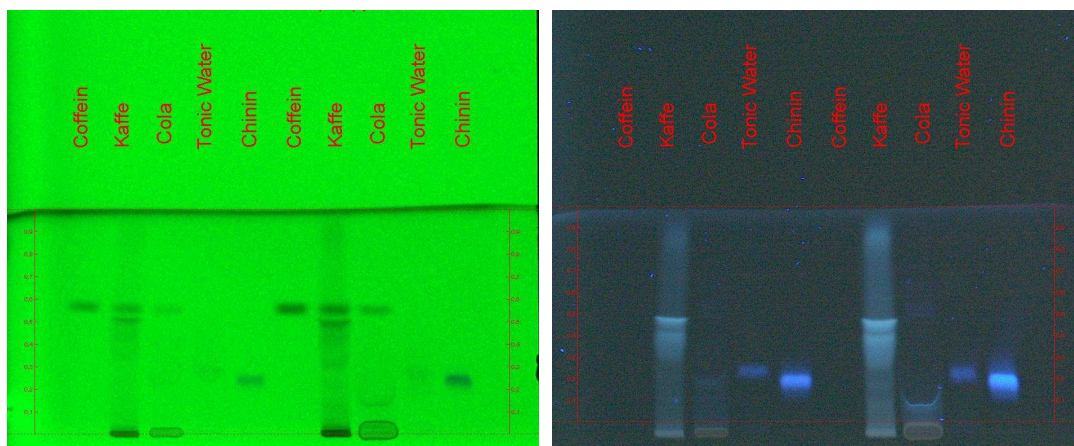




Detection of Caffeine in Coffee, Coke and Quinine in Tonic Water on TLC Silica gel 60 G F₂₅₄

Chromatographic Conditions

Plate: TLC Silica gel 60 G F₂₅₄ 10x10 cm
 Article number: 1.00390.0001
 Mobile Phase (v/v): Isopropanol / n-Heptan / Water 7:3:1 +1 % TEA
 Migration distance: 5 cm
 Chamber: normal chamber without chamber saturation
 Detection: UV 254 nm and 366 nm
 Sample preparation: Preparation of coffee in a coffee machine. 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	Caffeine	0.1	Methanol	1
2	Coffee extract			1
3	Coca Cola			1
4	Tonic water			1
5	Quinine	1	Methanol	1
6	Caffeine	0.1	Methanol	2
7	Coffee extract			2
8	Coca Cola			2
9	Tonic water			2
10	Quinine	1	Methanol	2