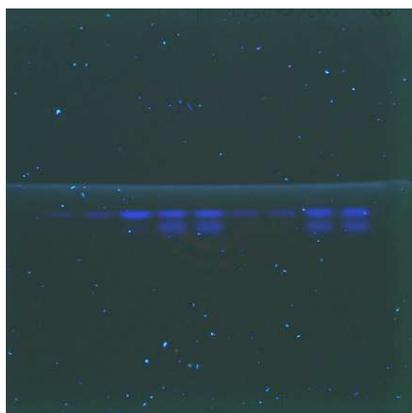




Determination of acetylsalicylic acid in painkillers on HPTLC Silica gel 60 F₂₅₄

Chromatographic Conditions

Plate: HPTLC Silica gel 60 F₂₅₄ 10x10 cm
 Article number: 1.05628.0001
 Mobile Phase (v/v): Ethylacetate (1.00868) / Methanol (1.06007) 4 : 1
 Migration distance: 5 cm
 Chamber: Normal chamber without chamber saturation
 Detection: 366 nm
 Sample preparation: 1 Tablet was dissolved in 10 ml Water and filtered over a 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, 2, 3	Acetylsalicylic acid	1	Water	0.5, 1.0, 2.0	84
4	Product 1		Water	2	84
5	Product 1		Water	2	84
6	Product 2		Water	2	84
7	Product 2		Water	2	84
8	Product 3		Water	2	84
9	Product 3		Water	2	84