

Separation of sugars on Silica gel 60

Plate	HPTLC Silica gel 60 20x10cm (P/N 1. 05641, alt. P/N 1. 05631)
Mobile phase (v/v)	1. 2-propanol/chloroform/methanol/ammonia25%/water (45/45/20/4/6) 2. water/ACN (15/85) 3. water/ACN (15/85) 4. water/ACN (15/85) 5. water/ACN (15/85)
Migration distance	Each 7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. β -Sitosterol 2. Diosgenin 3. Lanosterol (all 0,2%)
Application volume	2 μ l
Detection	Blei(IV)-acetat-Dichlorfluorescein-Reagent prepare a mixture of 3,75ml Blei(IV)-acetat saturated in glacial acetic acid and 3,75ml (0,22g Dichlorfluorescein in 100ml ethanol) and dilute to 150ml with toluene, spray 20s and heat 30min at 100°C, Detection UV @ 366nm

Chromatographic Data

No.	Compounds	Concentration (mg/ml)	Solvent
1	Raffinose	0,5	water
3	Maltose-Monohydrat	0,5	water
5	Saccharose	0,5	water
7	Galactose	0,5	water
9	D(+)-Glucose	0,5	water
11	D(-)-Fructose	0,5	water
13	D(-)-Arabinose	0,5	water
15	D(-)-Ribose	0,5	water
2,4,6,8,10,12,14,16	sugar mixture	0,5	water

