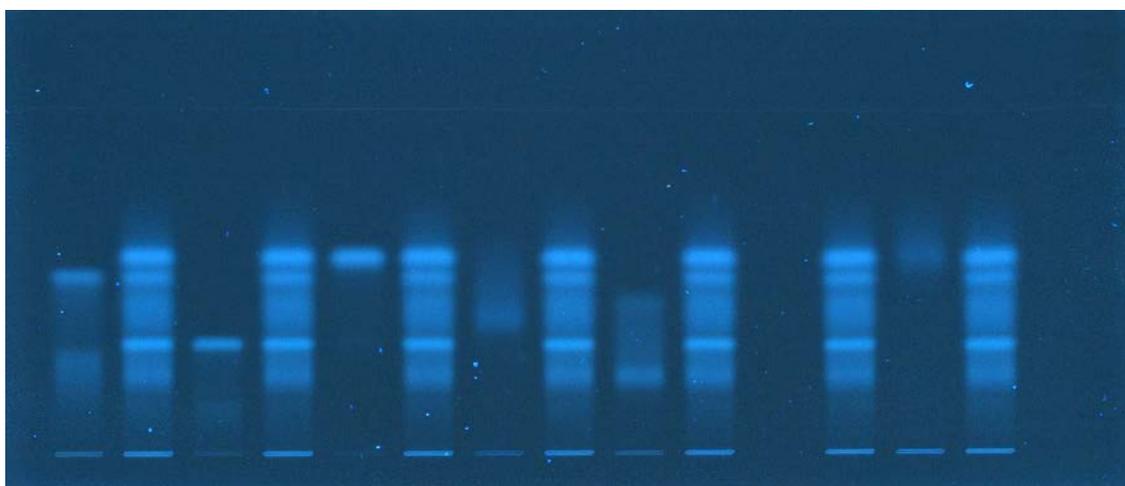


# Separation of sugars on Silica gel 60 NH<sub>2</sub>

Plate	HPTLC Silica gel 60 NH <sub>2</sub> 20x10cm (P/N 1. 13192)
Mobile phase (v/v)	1) water/ACN (30/70) 2) ethylacetate/pyridine/water/glacial acetic
Migration distance	1) 55mm, 2) 75mm
Application volume	2 µL
Chamber	Normal chamber without chamber saturation
Detection	Heat 5min at 150°C, Detection



## Chromatographic data

No.	Compounds	Concentration (mg/ml)	Solvent	hRf
1	D(+)-glucose	0,5	water	-
3	maltose-monohydrate	0,5	water	-
5	D(-)-fructose	0,5	water	-
7	D(-)-arabinose	0,5	water	-
9	D(+)-galactose	0,5	water	-
11	saccharose	0,5	water	-
13	D(-)-ribose	0,5	water	-
15	raffinose	0,5	water	-
2,4,6,8,10,12,14	sugar mixture	0,5	water	-