

# Separation of digested proteins on Silica gel 60 RP-8

Plate	HPTLC Silica gel 60 RP-8 F254s (P/N 1. 13725)
Mobile phase (v/v)	methanol/water (70/30) + 0,1mol/l ammonium
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Detection	0,02% fluorescamin solution in acetone spray 20s and dry 10min at room temperature Detection UV @ 366nm

## Chromatographic data

No.	Compounds	Concentration(mg/ml)	Solvent	Application volume(μl)
1	lysozyme digested	2,0	water/ACN 85/15	7,0
2	myoglobin digested	2,0	water/ACN 85/15	7,0
3	cytochrome c digested	2,0	water/ACN 85/15	7,0
4	β-casein digested	2,0	water/ACN 85/15	7,0
5	BSA digested	2,0	water/ACN 85/15	7,0

