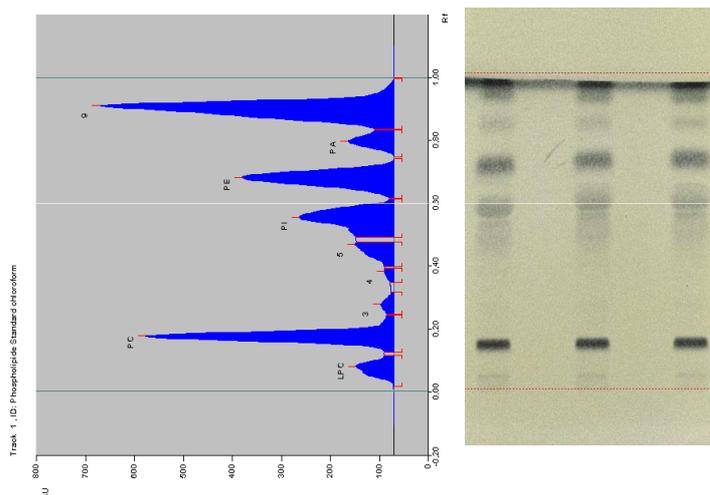




# Separation of Phospholipids on HPTLC Silica gel 60 F<sub>254</sub>

## Chromatographic Conditions

Phase: HPTLC Silica gel 60 F<sub>254</sub> 20x10 cm  
 Article number: 1.05642.0001  
 Mobile Phase (v/v): Chloroform / Methanol / Aceton / Water 18:15:3:1  
 Migration distance: 5 cm  
 Chamber: Normal chamber without chamber saturation  
 Derivatisation: Molybdätosphoric acid (1.00480.0100)  
 Detection: White light  
 Scan wave length: 720 nm



## Chromatographic Data

Track No.	Compounds	Concentration (mg/ml)	Solvent	Application volume $\mu$ l	hRF
1-3	Standard	1.0	Chloroform	5.0	
	Standard:				
	Triglyceride (TG)				95
	Fatty acid (FFA)				89
	Phosphatidic acid (PA)				73
	Phosphatidylethanolamine (PE)				61
	Phosphatidylinositol (PI)				54
	Phosphatidylcholine (PC)				14
	Lysophosphatidylcholine (LPC)				4