

Separation of nitroaniline, which have relative concentrations

HPTLC Si 60 F254

Plate	TLC aluminium sheets silica gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)		
Eluent	Carbon tetrachloride/acetic acid glacial 90/10 (v/v)		
Migration distance	6 cm		
Chamber	Normal chamber without chamber saturation		
Compounds	a.	b.	c.
1 p-Nitroaniline	1 %	1 %	1 %
2 m-Nitroaniline	0,005%	0,002%	0,001%
3 o-Nitroaniline	0,005%	0,002%	0,001%
Application volume	5 µl		
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm		

