

Separation of nitroaniline

HPTLC Si 60 F254

Plate	HPTLC pre-coated plate 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	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. p-Nitroaniline 2. m-Nitroaniline 3. o-Nitroaniline
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

