

Separation of phenylenediamines

HPTLC Si 60 F254

Plate	TLC pre-coated plate gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Chloroform/methanol, 90/10 (v/v) 80% saturated with water
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 1,4-Phenylenediamine 2. 1,3-Phenylenediamine 3. 1,2-Phenylenediamine (all 0,05%)
Application volume	1 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 280 nm

