

Separation of some chlorophenols

Plate	HPTLC pre-coated plate RP-2 F 254s (1.05747)
Eluent	Methanol/water 70/30 (v/v) with addition of 0,1 mol/l tetraethylammonium bromide
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Pentachlorophenol 2. 2,4,6-Trichlorophenol 3. 2,4-Dichlorophenol 4. 4-Chlorophenol
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

