

Separation of some chlorophenols

TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N 1. 15389, alt. P/N 1. 15423)
Eluent	Acetonitrile/water 70/30 (v/v)
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

