

# Separation of some chlorophenols

## TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N 1. 15388, alt. P/N 1. 15424)
Eluent	Methanol/water 65/35 (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

