

Separation of chlorophenols

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Methanol/water 90/10 (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 (all 0,1%)
Application volume	2 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

