

# Adsorption chromatographic separation of some chlorophenols

## HPTLC Silica gel 60 F254

Plate	HPTLC pre-coated plate silica gel 60 F 254 (P/N: 1.05642, Alternative P/N: 1.05548)
Eluent	Chloroform
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 2,4,6-Trichlorophenole (0,8%) 2. 2,4-Dichlorophenole (1,0%) 3. 4-Chlorophenole (1,0%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 224 nm

