

# Separation of chlorinated Hydrocarbons

## HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N: 1.05636, Alternative P/N: 1.12668)
Eluent	Petroleum benzine (40-60°)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Dieldrin (2%) 2. p,p'-DDT (2%) 3. Aldrin (3%)
Application volume	500 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 220 nm

