

Separation of some pesticides

HPTLC DIOL F254s

Plate	HPTIC pre-coated plate DIOL F 254s (P/N: 1.05636, Alternative P/N: 1.12668)
Eluent	Petroleum benzine (40 – 60°)/acetone 90/10 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Methoxuron 0,2% 2. Monuron 0,2% 3. Simazin 0,2% 4. Propazin 0,2% 5. Parathion-ethyl 0,2% 6. Alachlor 0,4% 7. Pendimethalin 0,2% 8. Trifluralin 0,2%
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

