

Separation of some pesticides

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Petroleum benzine (40–60°)/acetone 85/15 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Monuron 0,2% 2. Chlortoluron 0,2% 3. Simazin 0,2% 4. Sebuthylazin 0,2% 5. Propazin 0,2% 6. Metolachlor 0,4% 7. Pendimethalin 0,2% 8. Trifluralin 0,2%
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 230 nm.

