

Separation of pesticides

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Petroleum benzene (40–60°)/acetone 80/20 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	Derivatives of urea: 1. Mehoxuron 2. Monuron 3. Chlortoluron 4. Methobromuron (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm.

