

Separation of herbicides/fungicides

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 85/15 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Clopyralid (0,2%) 2. Metazachlor (0,4%) 3. Vinclozolin (0,4%) 4. Dinoseb(-acetate) (0,2%)
Application volume	500 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

