

# Separation of herbicides/fungicides

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F254s (1. 13192)
Eluent	Petroleum benzene (40-60°)/acetone 85/15 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Clopyralid 0,5% 2. Metazachlor 1,0% 3. Vinclozolin 0,4% 4. Dinoseb-(acetate) 0,2%
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

