

# Separation of herbicides/fungicides

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	n-hexane/toluene/acetone 70/26/4 (v/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

