

Separation of compounds of Aniline

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	n-hexane/toluene/acetone 70/25/5 (v/v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Metolachlor (1,0%) 2. Alachlor (2,0%) 3. Pendimethalin (0,2%) 4. Trifluralin (0,4%)
Application volume	100 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

