

Separation of derivatives of Urea

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Chloroform/toluene 80/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Methoxuron 2. Chlortoluron 3. Methobromuron (all 0,2%)
Application volume	150 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

