

Separation of derivatives of Urea

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N: 1.05636, Alternative P/N: 1.12668)
Eluent	Petroleum benzine (40–60°)/acetone 80/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Methoxuron 2. Monuron 3. Methobromuron (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 238 nm

