

# Separation of Phenoxy-carboxylic acids

## 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 85/15 (v/v)
Migration Distance	7 cm (2 x)
Chamber	Normal chamber without chamber saturation
Compounds	1. 2,4 - D 2. MCPA 3. Fenoprop (all 1%)
Application volume	750 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 220 nm

