

Separation of plasticizers

HPTLC Si 60 F254

Plate	TLC pre-coated plate gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Petroleum benzine 40–60°/acetone 90/10 (v/v)
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Dimethyl phthalate 2. Diethyl phthalate 3. Dibutyl phthalate (all 0,2%)
Application volume	1 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

