

# Separation of plasticizers

## HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Acetonitrile/methanol/water 45/45/10 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber with chamber saturation
Compounds	1. Dibutyl phthalate 2. Diethyl phthalate 3. Dimethyl phthalate 4. Phthalic acid (all 0,2%)
Application volume	1 $\mu$ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

