

# Separation of aromatic acids by ion-pair chromatography

## TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N: 1.15389, Alternative P/N: 1.15423)
Eluent	Acetone/water 50/50 (v/v) with addition of 0,1 mol/l TEABr
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 4-Methyl benzoic acid 2. Benzoic acid 3. 4-Hydroxy benzoic acid 4. 4-Amino benzoic acid 5. Naphthaline sulfonic acid (all 0,2%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

