

Separation of aromatic acids

TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N 1. 15388, alt. P/N 1. 15424)
Eluent	Acetonitrile/phosphate buffer 0,01 mol/l pH3 25/75 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 2-Methyl benzoic acid 2. Benzoic acid 3. Salicylic acid 4. 2-Nitro benzoic acid
Application volume	200 nl
Detection	In situ-evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

