

Separation of substituted benzoic acids by ion-pair chromatography

TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N: 1.15388, Alternative P/N: 1.15424)
Eluent	Acetone/water 40/60 (v/v) a. Without IPR b. With addition of 0,01 mol/l tetrapropylammonium bromide
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Benzoic acid 2. 2-Chloro benzoic acid 3. Salicylic acid
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

