

Separation of 4-Hydroxybenzoic acid esters

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	Aceton/water 40/60 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 4-Hydroxybenzoic acid propyl ester 2. 4-Hydroxybenzoic acid ethyl ester 3. 4-Hydroxybenzoic acid methyl ester 4. 4-Hydroxybenzoic acid
Application volume	300 nl
Detection	In situ-evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

