

The effect of stabilizer in the eluent on the separation of 4-Hydrobenzoic acid and some of its esters

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s (1. 13724)
Eluent	Cyclohexane/chloroform 30/70 (v/v) a. Chloroform stabilized with ethanol b. Chloroform stabilized with 2-methyl-2-buten
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. PHB 2. PBH-methylester 3. PHB-ethylester 4. PHB-propylester (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

