

Separation of preservatives

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Ethanol/water/acetone 20/8g/10 (v/v/v) with addition of 0,1 mol/l tetraethylammonium chloride
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Benzoic acid 2,0% 2. Sorbic acid 0,1% 3. p-Hydroxybenzoic acid 0,1% 4. Gallic acid 0,1%
Application volume	100 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

