

Separation of benzoic acid and sorbic acid

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Ethanol/water 20/80 (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 (0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

