

# Separation of benzoic acid and sorbic acid

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Petroleum benzine 40-60°/acetone 65/35 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Digitoxigenin 2. Digoxigenin 3. Digoxin (all 0,05%)
Application volume	200 nl
Detection	Spray reagent: MnCl <sub>2</sub> - sulfuric acid heating up to 110° C for 5 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

