

Separation of phenols

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Ethanol/water 40/60 (v/v) with addition of 0,2 mol/l lithium chloride
Migration Distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. p-Nitrophenol (0,5%) 2. p-Cresol (2,0%) 3. Pyrocatechol (0,5%) 4. Hydroquinone (0,4%) 5. o-Toluidine (0,5%)
Application volume	500 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

