

Separation of biphenylols

HPTLC Si 60 F254

Plate	HPTLC pre-coated plate silica gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Toluene/chloroform/methanol 80/10/10 (v/v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 1,1'-Biphenyl-4,4'-diol 2. 1,1'-Biphenyl-2,2'-diol 3. 1,1'-Biphenyl-4-ol 4. 1,1'-Biphenyl-2-ol
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

