

Separation of biphenylols

TLC Si 60 F254

Plate	TLC pre-coated plate gel 60 F254 with concentrating zone (P/N 1. 11798, alt P/N 1. 05583)
Eluent	Toluene/chloroform/methanol 80/10/10 (v/v/v)
Migration distance	10 cm
Chamber	Normal chamber with chamber saturation
Compounds	1. 1,1'-Biphenyl-4,4'-diol 0,01% 2. 1,1'-Biphenyl-2,2'-diol 0,05% 3. 1,1'-Biphenyl-4-ol 0,01% 4. 1,1'-Biphenyl-2-ol 0,05% 5. Biphenyl 0,1%
Application volume	5 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

