

Separation of some polyvalent phenols

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Chloroform/methanol 50/50 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Resorcinol 2. Hydroquinone 3. l-Naphthol 4. m-Cresol
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

