

Separation of some polyvalent phenols

HPTLC NH₂F 254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Acetone/chloroform 50/50 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Phloroglucinol 2. Resorcinol 3. Hydroquinone 4. 1-Naphthol 5. m-Cresol (all 0,2%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

