

Separation of some benzenecarboxylic acids

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N 1. 13192, alt. P/N 1. 15647)
Eluent	Ethanol/ammonia pH 12 60/40 (v/v) with the addition of 0,18 mol/l NaCl
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Pyromellitic acid 2. Benzene-1,2,3-tricarboxylic acid 3. Benzene-1,2,4-tricarboxylic acid 4. Phthalic acid 5. Benzoic acid (all 0,2%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm.

