

Separation of some naphthalene sulfonic acids

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Ethanol/ammonia pH 12 60/40 (v/v) with addition of 0,18 mol/l NaCl
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Naphthalene-1,3,7-trisulfonic acid 2. Naphthalene-1,3,6-trisulfonic acid 3. Naphthalene-1,5-disulfonic acid 4. Naphthalene-1-sulfonic acid (all 0,5%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

