

Separation of oligo nucleotides

HPTLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (1.13724)
Eluent	Methanol/buffer solution pH3 70/30 (v/v) with addition of 0.1 mol/l tetramethylammonium chloride
Migration Distance	12 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. ApApA 2. ApApC 3. ApApU 4. ApUpG (all 0,1%)
Application volume	500 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

