

Separation of oligo nucleotides

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254 (1.13724)
Eluent	Methanol/water 60/40 (v/v) with addition of 0,1 mol/l tetramethylammoniumbromide
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Benzoic acid 2. Isophthalic acid 3. Phthalic acid (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

