

Separation of oligo nucleotides

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Ethanol/water 10/90 (v/v) with addition of 0,1 mol/l lithium chloride
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. ApUpG 2. ApApU 3. ApApC (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

