

# Separation of oligo nucleotides

## HPTLC NH2 F254s

|                    |                                                                            |
|--------------------|----------------------------------------------------------------------------|
| Plate              | HPTLC pre-coated plate NH2 F 254s (P/N: 1.13192, Alternative P/N: 1.15647) |
| Eluent             | Ethanol/water 60/40 (v/v) with addition of 0,2 mol/l lithium chloride      |
| Migration Distance | 7 cm                                                                       |
| Chamber            | Normal chamber without chamber saturation                                  |
| Compounds          | 1. ApUpG<br>2. ApApU<br>3. ApApC<br>4. ApApA (all 0,1%)                    |
| Application volume | 300 nl                                                                     |
| Detection          | In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm                |

