

Separation of N-compounds on TLC pre-coated plate RP-8 F 254s by ion-pair chromatography

TLC Silica gel 60 RP-8 F254s

| | |
|--------------------|--|
| Plate | TLC pre-coated plate RP-8 F 254s (P/N: 1.15388, Alternative P/N: 1.15424) |
| Eluent | Acetone/water 40/60 (v/v) with addition of 0,1 mol/l octane-1 sodium sulfonate |
| Migration Distance | 7 cm |
| Chamber | Normal chamber without chamber saturation |
| Compounds | 1. Eupaverine chloride 2. Papaverine 3. Codeine 4. Caffeine |
| Detection | In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm |

