

Separation of alkaloids by ion-pair chromatography

TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N: 1.15389, Alternative P/N: 1.15423)
Eluent	Methanol/aqueous solution of 0,05 mol/l KH_2SO_4 pH4 and 0,1 mol/l octane-1-sodium sulfonate 40/60 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Eupaverine chloride 2. Papaverine 3. Caffeine 4. Codeine
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

