

Separation of morphine derivatives by ion-pair chromatography

TLC Silica gel RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N: 1.15389, Alternative P/N: 1.15423)
Eluent	Acetone/water 40/60 (v/v) with addition of 0,1 mol/l heptanesulfonic acid sodium salt
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Dionin (Ethylmorphine) 2. Codeine (Methylmorphine) 3. Morphine hydrochloride
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

