

Separation of tropane-alkaloids

TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (P/N 1. 15389, alt. P/N 1. 15423)
Eluent	Aceton/water 40/60 (v/v) with addition of 0.3 mol/l NaCl
Migration distance	7 cm
Chamber	Normal chamber with chamber saturation
Compounds	1. Apotatropine 2. Atropine 3. Scopolamine
Detection	In situ-evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

