

Separation of alkaloids

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Methanol/l-propanol /water/ammonia solution 25% 30/20/50/1 with addition of 0,1 mol/l ammonium bromide
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Emetine chloride (0,4%) 2. Quinine (0,1%) 3. Atropine (3,0%) 4. Brucine (0,1%) 5. Scopolamine (3,0%) 6. Caffeine (0,3%) 7. Nicotinamide (0,2%) 8. DL-Tropic acid (3,0%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

