

# Separation of some drugs

## TLC Silica gel 60 RP-18 F254s

Plate	TLC pre-coated plate RP-18 F 254s (E.Merck) (P/N: 1.15389, Alternative P/N: 1.15423)
Eluent	Acetone/water 50/50 (v/v) with addition of 0,1 mol/l tetraethylammonium bromide
Chamber	Normal chamber without chamber saturation
Compounds	1. Methaqualone 2. Phenobarbital 3. Diphenhydramine 4. Barbitol 5. Codeine 6. Morphine
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

