

Separation of some cholesteryl esters

HPTLC Si 60 F254

Plate	HPTLC pre-coated plate silica gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Toluene/petroleum benzene 40-60° 40/60 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Cholesteryl butonate 2. Cholesteryl hexonate 3. Cholesteryl linoleate 4. Cholesteryl oleate 5. Cholesteryl erucate (all 0,01%)
Application volume	5 µl
Detection	Spray reagent: MnCl ₂ -sulfuric acid heating up to 110°C for 5 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

