

Separation of cholesterol and its bile acid metabolites

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	Methanol/acetone/acetic acid 85/10/5 (v/v/v)
Migration distance	10 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Cholesterol 2. 7-Hydroxycholesterol 3. Lithocholic acid 4. Cholic acid methyl ester 5. Cholic acid 6. Dehydrocholic acid
Application volume	400 nl
Detection	Spray reagent: MnCl_2 – sulfuric acid heating up to 120°C for 120 min. UV 366 nm.

