

Separation of fatty acids as isatinmethyl esters

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s (1. 13724)
Derivatisation	Reaction of the fatty acids with 1-chloromethyl-isatin
Eluent	Methanol/water 88/12 (v/v)
Chamber	Normal chamber without chamber saturation
Compounds	Isatin esters of: 1. Myristic acid 2. Tridecanoic acid 3. Lauric acid 4. n-Capric acid
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 313 nm
Application of reaction mixture	a. Chromatogram of reaction mixture without fatty acids b. Chromatogram of reaction mixture

