

Separation of fatty acids

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 254s F (1. 13724)
Eluent	Ethanol/water 75/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Palmitic acid (8%) 2. Myristic acid (8%) 3. Lauric acid (8%) 4. Capric acid (8%) 5. Caprylic acid (8%)
Application volume	200 nl
Detection	Dipped in dichlorofluoresceince (0,2% in ethanol)

