

# Separation of edible fats

## HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s (1. 13724)
Eluent	Acetone/acetonitrile/Toluene/Chloroform/n-Hexane 50/35/5/5/5 v/v
Migration Distance $z_f$	7.5 cm
Chamber	Normal chamber without chamber saturation
Application volume	20 $\mu$ l (10mg/5ml)
Detection	Dipped in 1% molybdato-phosphoric acid in methanol, heating up to 110°C

