

# Separation of fats and oils

## HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Acetone/acetonitrile/chloroform 50/40/20 (v/v/v)
Migration Distance	7,5 cm
Chamber	Normal chamber with chamber saturation
Application volume	5 µl (10mg/5ml)
Detection	Dipped in 1% molybdatophosphoric acid in methanol heating up to 110°C

