

# Separation of cholesterol and sitosterol

Plate	HPTLC pre-coated plate RP-18 with concentrating zone (1.15498)
Eluent	Cyclohexane/ethyl acetate/acetonitrile 60/30/10 (v/v/v) Temperature of development: +10°C
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Cholesterol 2. Sitosterol (all 0,02%)
Application volume	5 µl
Detection	MnCl <sub>2</sub> - sulfuric acid (spray reagent) heating up to 120° C UV 366 nm

