

Separation of steroids

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	N-hexane/acetone 90/10 (v/v) temperature of development: + 4°C
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. β -Sitosterol 2. Diosgenin 3. Lanosterol (all 0,2%)
Application volume	100 nl
Detection	Spray reagent: MnCl_2 -sulfuric acid heating up to 110°C for 5 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

