

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/acetic acid glacial 90/10/1 (v/v/v) with addition of 0,1 mol/l lithium chloride
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Quercetin 2. Penogenin 3. Diosgenin 4. Stigmasterol 5. Lanosterol (all 0,1%)
Application volume	200 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.

