

Separation of testosterones

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Di-iso-propyl ether
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 1,4 α -Androstatrien-3, 17-dione 2. Epiandrosterone 3. 5 β -Dihydrotestosterone 4. Androstane-3, 17-dione (all 0,1%)
Application volume	200 nl
Detection	Spray reagent: MnCl ₂ -sulfuric acid heating up to 110°C for 10 min In-situ evaluation with TLC/HPTLC scanner (Camag) λ 540 nm

