

Separation of androgens

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Petroleum benzine (40-60°)/acetone 80/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Testosterone 2 5-Dehydro-Androsterone 3 Androsterone (all 0,2%)
Application volume	200 nl
Detection	Spray reagent: 5% methanolic perchloric acid heating up to 110°C for 5 min. UV 366 nm

