

Separation of steroids

HPTLC Silica gel 60 RP-8 F254s

Plate	HPTLC pre-coated plate RP-8 F 254s (1. 13725)
Test compounds	1. Methyltestosterone 2. Reichstein's substance S 3. Hydrocortisone
Application volume	200 nl (10mg/10ml of each compound)
Eluent	Acetone/water 60/40 v/v
Migration distance z_f	5 cm
Type of chamber	Normal chamber without chamber saturation
Detection	UV 254nm, TLC/HPTLC Scanner from Camag

