

# Separation of steroids

## HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s (1. 13724)
Eluent	Acetone/water 60/40 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Methyltestosterone 2. Reichstein's substance S 3. Hydrocortisone (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

