



Separation of hormones (drugs for the climacteric period)

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1.13124)
Eluent	Acetonitrile/water 60/40 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Testosterone 1,6% 2. Estradiol 0,06%
Application volume	5 µl
Detection	Dipped in MnCl ₂ -sulfuric acid heating up to 110° C for 2 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

