

# Separation of hormones (drug for the climacteric period)

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1. 13124)
Eluent	Acetonitrile/water 40/60 (v/v)
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Estradiol 0,4% 2. Estriol 0,2%
Application volume	5 $\mu$ l
Detection	Dipped in MnCl <sub>2</sub> -sulfuric acid heating up to 110°C for 2 min. In-situ evaluation with TLC/HPTLC scanner (Camag) wave-length 440 nm.

