

# Separation of estradiol and some of its esters

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1. 13124)
Eluent	Acetonitrile/water 75/25 (v/v)
Migration distance	6 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Estradiol 17 $\beta$ -cypionate 2. Estradiol 17 $\beta$ -valerate 3. Estradiol 17 $\beta$ -acetate 4. Estradiol (all 0,0125%)
Application volume	5 $\mu$ l
Detection	Dipped in MnCl <sub>2</sub> -sulfuric acid heating up to 110°C for 10 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm.

