

Separation of estrone derivatives

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Petroleum benzine 40–60°C/acetone 80/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Estriol 2. Estradiol 3. Estrone (all 0,1%)
Application volume	200 nl
Detection	Spray reagent: MnCl ₂ -sulfuric acid heating up to 120° C for 10 min. UV 366 nm

