

Separation of weak polar steroids

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Di-iso-propyl ether/acetone/acetic acid glacial 80/20/1 (v/v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Dexamethasone 2. Cortisol 3. Corticosterone 2. Deoxycorticosterone (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

