

Separation of anabolic compounds

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F 254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Di-iso-propyl ether/acetic acid glacial 100/1 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 19-Nortestosterone 2. Medroxyprogesterone 3. Progesterone 4. α -Dienestrol (all 0,1%)
Application volume	300 nl
Detection	Spray reagent: MnCl_2 – sulfuric acid heating up to 110°C for 5 min. In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

