

# Separation of progesterones

## HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Petroleum benzine (40-60°)/acetone 75/25 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Prednisone 2. 11,17-Ketoprogesterone 3. Reichstein's substance "S" 4. Progesterone 5. Pregnadienoloneacetate (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

