

# 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	<ol style="list-style-type: none"> <li>1. 9<math>\alpha</math>-Flourohydrocortisone</li> <li>2. Allo-Dihydrocortisone</li> <li>3. 11<math>\alpha</math>-Pregnane-3<math>\alpha</math>-2<math>\alpha</math>-diol</li> <li>4. Pregnanediol</li> <li>5. Pregnadienoloneacetate</li> <li>6. Progesterone</li> <li>7. Pregnadienoloneacetate (all 0,1%)</li> </ol>
Application volume	300 nl
Detection	<p>Spray reagent: MnCl<sub>2</sub>-sulfuric acid heating up to 110°C for 5 min.</p> <p>In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm.</p>

