

# 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. 16-Methylprednisolone</li> <li>2. 16-Methyl-4-pregnene-3,20-dione-11<math>\alpha</math>-17<math>\alpha</math>-diol</li> <li>3. 11<math>\alpha</math>-Hydroxyprogesterone</li> <li>4. Deoxycorticosterone</li> <li>5. 6-Methyl-6-dehydro-17<math>\alpha</math>-progesterone</li> <li>6. Progesterone</li> <li>7. Pregnadienoloneacetate (all 0,1%)</li> </ol>
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

