

Separation of digitalis glycosides

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate DIOL F 254s (P/N 1. 05636, alt. P/N 1. 12668)
Eluent	Ethyl acetate/ammonia solution 25% 100/1 (v/v)
Migration distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	<ol style="list-style-type: none"> 1. Lanatoside C 2. Digoxin 3. Digitoxin 4. Digoxigenin 5. α-Acetyldigoxin 6. Digitoxigenin (all 0,1%)
Application volume	100 nl
Detection	<p>Spray reagent: $MnCl_2$ – sulfuric acid heating up to 110°C for 10 min.</p> <p>In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm</p>

