

Separation of mycotoxins

TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N 1. 15388, alt. P/N 1. 15424)
Eluent	Acetonitrile/buffer solution pH 2,00 50/50 (v/v)
Migration distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Zearalenone 2. Zeralenol 3. Rubratoxin B 4. Penicillic acid 5. Patulin (all 0,1%)
Application volume	300 nl
Detection	In situ-evaluation with TLC/HPTLC scanner (Camag) UV 248 nm

