

Separation of mycotoxins

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Acetonitrile/buffer solution pH 2,00 50/50 (v/v)
Migration Distance	7 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	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

