

Separation of sulfonamides

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1.13124)
Eluent	Methanol/water 50/50 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Sulfamethoxazol 0,4% 2. Trimethoprim 0,08%
Application volume	5 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

