

Separation of sulfonamides

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Acetonitrile/methanol/water 20/20/60 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Trimethoprin 2 Sulfadimethoxine 3 Sulfamethoxazole 4 Sulfathiazole 5 Sulfanilamide (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

