

# Separation of sulfonamides

Plate	HPTLC pre-coated plate RP-18WF 254s with concentrating zone (1.13124)
Eluent	Acetone/water 50/50 (v/v) with addition of 0,1 mol/l TEACl
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Sulfadiazine 0,41% 2. Trimethoprim 0,09%
Application volume	5 $\mu$ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

