

# Separation of ingredients in analgesic agents

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1.13124)
Eluent	Acetonitrile/water 40/60 (v/v) with addition of 0,1 mol/l ammonium bromide
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Acetylsalicylic acid 0,325% 2. Paracetamol 0,250%
Application volume	5 $\mu$ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

