

# Separation of ingredients in analgesics

## HPTLC Silica gel 60 RP-18 WF254s

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Acetonitrile/water 40/60 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Phenacetin (0,2%) 2. Caffeine (0,3%) 3. Acetylsalicylic acid (0,5%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 248 nm

