

Separation of ingredients in analgesics

HPTLC Silica gel 60 RP-18 WF254s

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Methanol/buffer solution pH 7/water 60/20/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Codeine (0,35%) 2. Phenacetin (0,10%) 3. Acetylsalicylic acid (0,50%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

