

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)
Migration distance	6 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Propyphenazone 0,15% 2. Phenacetin 0,25% 3. Caffeine 0,03%
Application volume	5 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 248 nm.

