

Separation of ingredients in anti-asthmatic agents

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1. 13124)
Eluent	Acetonitrile/water 30/70 (v/v) with addition of 0,1 mol/l TEACl
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Theophylline 0,25% 2. Ephedrine 0,125%
Application volume	5 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 210 nm.

