

Separation of ingredients in circulatory stimulants

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1. 13124)
Eluent	Aceton/water 30/70 (v/v) with addition of 0,1 mol/l ammonium bromide
Migration distance	7 cm
Chamber	Normal chamber with chamber saturation
Compounds	1. Papaverine hydrochloride 0,003% 2. Paracetamol 0,1% 3. Theobromine 0,05%
Application volume	5 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 248 nm.

