

Separation of ingredients in dermatics

HPTLC Silica gel 60 RP-18 WF254s

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Chloroform/methanol 95/5 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber with chamber saturation
Compounds	1. Prednisolone 2. Linolic acid (all 0,2%)
Application volume	200 nl
Detection	Spray reagent: MnCl_2 - sulfuric acid heating up to 100°C for 10 min. In-situ evaluation with TLC/HPTLC scanner (Camag) λ 500 nm

