

Separation of vitamine

TLC aluminium oxide 150 F254

Plate	TLC pre-coated plate aluminium oxide 150 F 254 (type T) (1.05551)
Eluent	Chloroform/methanol 50/50 (v/v) 70% saturated with 1N acetic acid
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Vitamin B1 0,5% 2. Vitamin B12 0,1% 3. Nicotinamide 0,5%
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

