

Separation of water soluble vitamins

HPTLC Silica gel 60 RP-18 F254s

Plate	HPTLC pre-coated plate RP-18 F 254s with concentrating zone (1.15498)
Eluent	Acetonitrile/dichlormethane/water 90/10/10 (v/v) with addition of 0,1 mol/l lithium chloride
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Nicotin amide 2. Iso-Nicotinic acid 3. Nicotinic acid 4. L-Ascorbic acid (all 0,05%)
Application volume	1 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

