



# Separation of vitamine

## HPTLC NH<sub>2</sub> F254s

Plate	HPTLC pre-coated plate NH <sub>2</sub> F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Acetonitrile/water 70/30 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Vitamin B6 2. Vitamin B12 3. Vitamin B1
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 280 nm

