

# Separation of PTH amino acids

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Methanol/l-propanol /water 25/20/55 (v/v/v) with addition of 0,1 mol/l lithium chloride
Migration Distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. PTH-OL-Tryptophan 2. PTH-OL-Isoleucine 3. PTH-OL-Tyrosine 4. PTH-Glycine 5. PTH-OL-Arginine 6. PTH-OL-Cysteic acid (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

