

# Separation of PTH amino acids

## HPTLC CN F254s

|                    |                                                                                                                     |
|--------------------|---------------------------------------------------------------------------------------------------------------------|
| Plate              | HPTLC pre-coated plate CN F 254s (1.16464)                                                                          |
| Eluent             | Methanol/1-propanol/water 30/20/50 (v/v/v) with addition of 0,1 mol/l lithium chloride                              |
| Migration Distance | 7 cm                                                                                                                |
| Chamber            | Normal chamber without chamber saturation                                                                           |
| Compounds          | 1. PTH-DL-Lysine<br>2. PTH-DL-Valine<br>3. PTH-DL-Proline<br>4. PTH-DL-Glutamic acid<br>5. PTH-DL-Serine (all 0,2%) |
| Application volume | 200 nl                                                                                                              |
| Detection          | In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm                                                         |

