

# Separation of dansyl amino acids

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Methanol/l-propanol/water 50/20/30 (v/v/v) with addition of 0,1 mol/l lithium chloride
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Dansyl-L-tyrosine 2. Dansyl-L-lysine 3. Dansyl-L-phenylalanine 4. Dansyl-L-proline 5. Dansyl-L-serine (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 366 nm

