

Separation of DNP amino acids

HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F 254s (1.16464)
Eluent	Ethanol/acetone/water 25/30/45 (v/v/v) with addition of 0,2 mol/l lithium chloride
Migration Distance	8 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Ne-DNP-L-Lysine hydrochloride 2. N,N'-di-DNP-L-Cystine 3. N-DNP-DL-Methionine 4. N-DNP-L-Alanine 5. N-DNP-L-Glutamic acid cyclohexylamine salt 6. N-DNP-L-Asparagine (all 0,2%)
Application volume	400 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

