

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. N,O-DNP-L-Tyrosine 2. N,S-di-ONP-L-Cysteine 3. N-DNP-L-Isoleucine 4. N-DNP-Glycine 5. N-DNP-L-Serine 6. N-DNP-L-Arginine (all 0,2%)
Application volume	400 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

