

# Separation of Nitrophenols

## HPTLC CN F254s

Plate	HPTLC pre-coated plate CN F254s (1.16464)
Eluent	Carbon tetrachloride/methanol 48/2 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 2,3-Dinitrophenol 2. 4-Nitrophenol 3. 3-Nitrophenol 4. 2,6-Dinitrophenol 5. 2,4-Dinitrophenol 6. 2,5-Dinitrophenol 7. 2-Nitrophenol (all 0,2%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

