

# Separation of dinitrophenols

## HPTLC RP 18W F254s

|                    |   |
|--------------------|---|
| Plate              | HPTLC pre-coated plate RP 18W F254s (1.13724)   |
| Eluent             | Acetone/water 30/70 (v/v)   |
| Migration Distance | 7 cm  |
| Chamber            | Normal chamber without chamber saturation   |
| Compounds          | 1. 2,3-Dinitrophenol<br>2. 2,5-Dinitrophenol<br>3. 2,4-Dinitrophenol<br>4. 2,6-Dinitrophenol (all 0,2%) |
| Application volume | 100 nl  |
| Detection          | In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm   |

