

Separation of uridine and its phosphates

TLC PEI-Cellulose F

Plate	TLC pre-coated plate PEI-Cellulose F (P/N: 1.05725, Alternative P/N: 1.05579)
Eluent	Acetic acid 1N/ethanol 80/20 (v/v) with addition of 0,5 mol/l lithium chloride
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. UTP 2. UDP 3. UDP-Glucose 4. UMP 5. Uridine (all 0,1%)
Application volume	400 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag)UV 270 nm

