

Separation of nucleosides

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1.13124)
Eluent	Water/ethanol 95/5 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Adenosine 2. Guanosine 3. Inosine 4. Cytidine 5. Uridine (all 0,005%)
Application volume	5 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

