

Separation of nucleosidemonophosphates

Plate	HPTLC pre-coated plate RP-18 WF 254s with concentrating zone (1.13124)
Eluent	Water with addition of 0,2 mol/l l-butanefulfonic acid sodium salt
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. AMP 2. GMP 3. IMP (all 0,01%)
Application volume	5 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 260 nm

