

Separation of nucleobases

Plate	HPTLC pre-coated plate RP-2 F 254s (1. 05747)
Eluent	Acetonitrile/water 60/40 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Cytosine 2. Adenine 3. Guanine 4. Uracil (all 0,1%)
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

