

Separation of nucleobases

HPTLC Silica gel 60 RP-18 WF254s

Plate	HPTLC pre-coated plate RP-18w F 254s (1.13124)
Eluent	Acetonitrile/water 60/40 (v/v)
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Cytosine 2. Adenine 3. Guanine 4. Uracil (all 0,1%)
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

