

Separation of nucleobases

HPTLC Silica gel 60 RP-2 F254s

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
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 264 nm

