

Separation of some pyrimidine bases

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Chloroform/methanol 60/40 (v/v) 80% saturated with water
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Uracil 2. Cytosine 3. Thymine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

