

Separation of some pyrimidine derivatives

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Chloroform/methanol 50/50 (v/v) 80% saturated with water
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. 2-Thiouracil 2. Uracil 3. Cytosine 4. Thymine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

