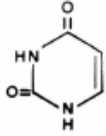
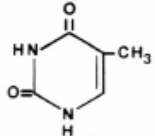
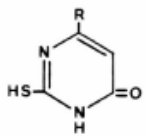


Separation of pyrimidine derivatives

HPTLC Si 60 F254

Plate	TLC pre-coated plate gel 60 F254 with concentrating zone (P/N 1. 13728, alt P/N 1. 13727)
Eluent	Chloroform/methanol 90/10 (v/v)
Migration distance	5,5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Uracil 2. Thymine 3. Thiouracil 4. 4-Methyl-2-thiouracil 5. 4-Propyl-2-thiouracil (All 0,01%)    R = -H R = -CH₃ R = -C₃H₇
Application volume	2 µl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

