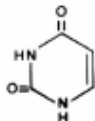
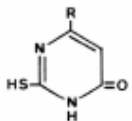


# Separation of pyrimidine derivatives

## 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 cm                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Chamber            | Normal chamber without chamber saturation                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                    |
| Compounds          | <ol style="list-style-type: none"> <li>1. Uracil</li> <li>2. Thiouracil</li> <li>3. 4-Methyl-2-thiouracil</li> <li>4. 4-Propyl-2-thiouracil</li> </ol> <p>(All 0,01%)</p> <div style="display: flex; align-items: center;"> <div style="margin-right: 20px;">  </div> <div>  <div style="margin-left: 10px;"> <p>R = - H</p> <p>R = - CH<sub>3</sub></p> <p>R = - C<sub>3</sub>H<sub>7</sub></p> </div> </div> </div> |
| Application volume | 2 µl                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                         |
| Detection          | In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                                  |

