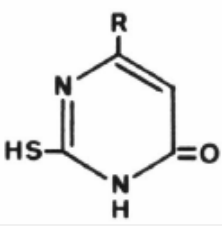


Separation of thiouraciles

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/2-propanol 90/10 (v/v)
Migration distance	6 cm
Chamber	Normal chamber without chamber saturation
Compounds	<p>All 0,01%</p> <div style="display: flex; align-items: center;"> <div style="margin-right: 20px;"> <p>Thiouracil $R = -H$</p> <p>4-Methyl-2-thiouracil $R = -CH_3$</p> <p>4-Propyl-2-thiouracil $R = -C_3H_7$</p> </div> <div>  </div> </div>
Application volume	2 μ l
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 265 nm

