

Separation of some purine derivatives

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Ethanol/water 80/20 (v/v) saturated with NaCl
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Uric acid 2. Xanthine 3. Hypoxanthine 4. Guanine 5. Adenine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 270 nm

