

Separation of xanthine and its N-methyl derivatives

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Methanol
Migration Distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Xanthine 0,03% 2. Theophylline 0,03% 3. Theobromine 0,06% 4. Caffeine 0,02%
Application volume	100 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

