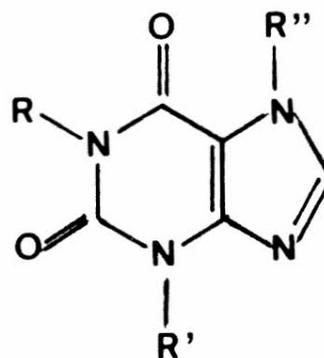
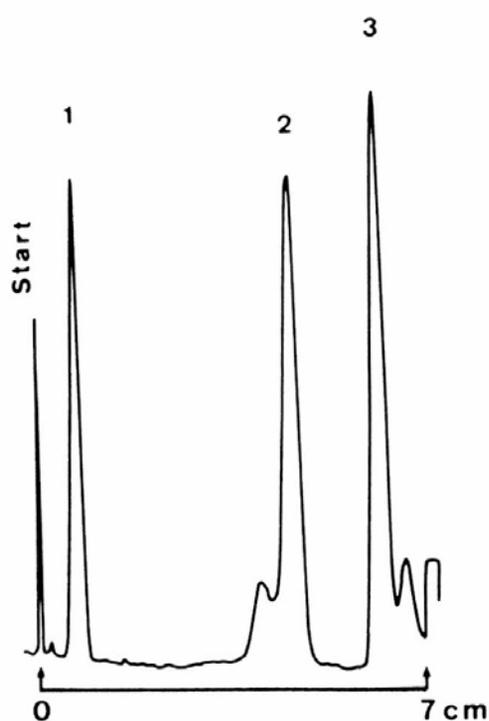


# Separation of the N-methyl derivatives of xanthine

## HPTLC NH<sub>2</sub> F254s

Plate	HPTLC pre-coated plate NH <sub>2</sub> F 254s (P/N: 1.13192, Alternative P/N: 1.15647)
Eluent	Methanol/acetonitrile/chloroform 10/80/10 (v/v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Theophylline 0,03% 2. Theobromine 0,06% 3. Caffeine 0,02%
Application volume	200 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm



R	R'	R''
CH <sub>3</sub>	CH <sub>3</sub>	H
H	CH <sub>3</sub>	CH <sub>3</sub>
CH <sub>3</sub>	CH <sub>3</sub>	CH <sub>3</sub>