

Separation of adenine derivatives

TLC Silica gel 60 RP-8 F254s

Plate	TLC pre-coated plate RP-8 F 254s (P/N 1. 15388 alt. P/N 1. 15424)
Eluent	Acetonitrile/water 20/80 (v/v)
Migration distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Adenine 2. Adenosine 3. cAMP 4. AMP (all 0,1%)
Application volume	300 nl
Detection	In situ-evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

