

Separation of adenine, adenosine and its phosphates

HPTLC DIOL F254s

Plate	HPTLC pre-coated plate NH ₂ F 254s (P/N 1. 13192, alt. P/N 1. 15647)
Eluent	Ethanol/water 30/70 (v/v) with addition of 0,2 mol/l NaCl
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. ATP 2. ADP 3. AMP 4. Adenine 5. Adenosine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm.

