

Separation of cytosine, cytidine and its phosphates

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Aceton/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. CTP 2. CDP 3. CMP 4. Cytosine, cytidine (all 0,1%)
Application volume	300 nl
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

