

Separation of guanine, guanosine and its phosphates

HPTLC NH₂ F254s

Plate	HPTLC pre-coated plate NH ₂ F254s (P/N 1. 13192, alt P/N 1. 15647)
Eluent	Methanol/water 5/95 (v/v) with addition of 0,2 mol/l NaCl
Migration distance	5 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. GTP 2. GDP 3. GMP 4. Guanine 5. Guanosine (all 0,1%)
Detection	In-situ evaluation with TLC/HPTLC scanner (Camag) UV 254 nm

