

Separation of phosphates

TLC Cellulose

Plate	TLC pre-coated plate cellulose (P/N: 1.05716, Alternative P/N: 1.05730)
Eluent	Dioxane / solution of 160g trichloroacetic acid, 8ml ammonia 25% per 1l in water 70/30 (v/v)
Migration Distance	10 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. $(\text{NaPO}_3)_3$ (1 %) 2. $\text{Na}_5\text{P}_3\text{O}_{10}$ (1 %) 3. $\text{Na}_4\text{P}_2\text{O}_7$ (1 %) 4. Na_2HPO_4 (0,2%)
Application volume	500 nl
Detection	Spray reagents (2xa) and then (1xb) a) $\text{Na}_2\text{MoO}_4 \cdot 2\text{H}_2\text{O}$ 18g NH_4NO_3 22,5g in 460ml water plus 40ml conc. HN03 b) $\text{Na}_2\text{S}_2\text{O}_5$ 150g Na_2SO_3 5g p-methylaminophenol 1g in 500ml water

