

Separation of amino acids

HPTLC Cellulose

Plate	HPTLC pre-coated plate cellulose without fluorescent indicator (P/N: 1.05651, Alternative P/N: 1.16092)
Eluent	l-propanol/water 80/20 (v/v)
Migration Distance	7 cm
Chamber	Normal chamber without chamber saturation
Compounds	1. Arginine 2. β -Alanine 3. α -Alanine 4. DL-Valine 5. L-Leucine
Detection	Spray reagent: ninhydrin heating up to 110°C for 5 min

