


Certificate of Analysis

Product Name:	Select-HA HiLadder				
Product Number:	HYA-HILAD				
Lot Number:	XCM090924				
Product Description:	Lyophilized white solids containing 5 Select-HA markers with molecular weight range between approximately 500 kDa and 1500 kDa				
Quality Assay:	<ul style="list-style-type: none"> • Electrophoresis • Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC) 				
Quality Parameters:	Parameter	Actual	Required	Pass	
	A. Markers Clarity on Gel	Clear	Clear	✓	
	B. Polydispersity				
	<u>Marker</u>	<u>Mn</u>	<u>Mw</u>	<u>Mw/Mn</u>	<u>Required</u> <u>Pass</u>
	#1	1271 kDa	1272 kDa	1.001	≤1.05 ✓
	#2	932 kDa	932 kDa	1.000	≤1.05 ✓
	#3	762 kDa	763 kDa	1.001	≤1.05 ✓
	#4	547 kDa	547 kDa	1.000	≤1.05 ✓
	#5	468 kDa	480 kDa	1.026	≤1.05 ✓
	Figure 1. Electrophoresis Analysis 5 µL of reconstituted Select-HA HiLadder (Lot# XCM090924; 1 vial reconstituted in 100 µL water) was run on 0.7% (w/v) agarose gel and stained with 0.005% Stains-all in 50% ethanol.				Marker #1 – 1272 kDa Marker #2 – 932 kDa Marker #3 – 763 kDa Marker #4 – 547 kDa Marker #5 – 480 kDa

This product is for research use only. Not intended or approved for diagnostic or therapeutic use.

Passes All Functional Specifications:  Angie M. Branch, Biology Manager

Date: 8/16/2024