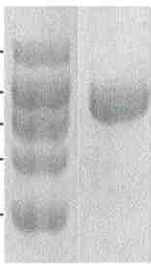
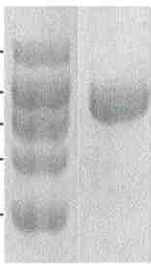
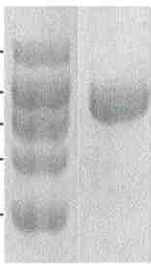


Echelon Biosciences Inc.

Certificate of Analysis

Product Name:	1000 kDa Select-HA																																						
Product Number:	HYA-1000																																						
Lot Number:	FVC070722B																																						
Product Description:	Lyophilized white solids containing 1 mg Select-HA (determined by Carbazole Assay)																																						
Quality Assay:	<ul style="list-style-type: none">• Electrophoresis• Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC)• Chromogenic Limulus Amebocyte Lysate (LAL) Assay																																						
Quality Parameters	<table border="1"><thead><tr><th>Parameter</th><th>Required</th><th>Actual</th><th>Pass</th></tr></thead><tbody><tr><td>A. Clarity on Gel</td><td>Clear</td><td>Clear</td><td>✓</td></tr><tr><td>B. Polydispersity</td><td>≤1.05</td><td>1.001</td><td>✓</td></tr><tr><td>C. Molecular Mass</td><td>950-1100 kDa</td><td>1062 kDa</td><td>✓</td></tr><tr><td>D. Endotoxin Level</td><td>≤0.1 EU/mg</td><td>≤0.1 EU/mg</td><td>✓</td></tr></tbody></table> <table><tr><td></td><td>1</td><td>2</td></tr><tr><td>1650 kDa —</td><td></td><td></td></tr><tr><td>1140 kDa —</td><td></td><td></td></tr><tr><td>918 kDa —</td><td></td><td></td></tr><tr><td>667 kDa —</td><td></td><td></td></tr><tr><td>509 kDa —</td><td></td><td></td></tr></table> <p>Electrophoresis Analysis Select-HA was run on 1% (w/v) agarose gel and stained with 0.005% Stains-All in 50% ethanol.</p> <p>Lane 1 – Select-HA HiLadder (HYA-HILAD-20, lot# HL201812)</p> <p>Lane 2 – 1000 kDa Select-HA (HYA-1000, lot# FVC070722B)</p>	Parameter	Required	Actual	Pass	A. Clarity on Gel	Clear	Clear	✓	B. Polydispersity	≤1.05	1.001	✓	C. Molecular Mass	950-1100 kDa	1062 kDa	✓	D. Endotoxin Level	≤0.1 EU/mg	≤0.1 EU/mg	✓		1	2	1650 kDa —			1140 kDa —			918 kDa —			667 kDa —			509 kDa —		
Parameter	Required	Actual	Pass																																				
A. Clarity on Gel	Clear	Clear	✓																																				
B. Polydispersity	≤1.05	1.001	✓																																				
C. Molecular Mass	950-1100 kDa	1062 kDa	✓																																				
D. Endotoxin Level	≤0.1 EU/mg	≤0.1 EU/mg	✓																																				
	1	2																																					
1650 kDa —																																							
1140 kDa —																																							
918 kDa —																																							
667 kDa —																																							
509 kDa —																																							

Passes All Functional Specifications:



Date: 8/3/22

Angie M. Branch, Biology Manager

For research use only. Not intended or approved for diagnostic or therapeutic use.