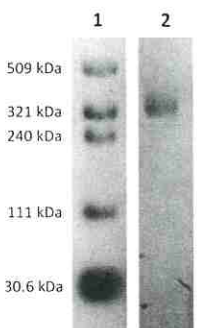


# Echelon Biosciences Inc.

## Certificate of Analysis

<b>Product Name:</b>	400 kDa Select-HA																						
<b>Product Number:</b>	HYA-0400																						
<b>Lot Number:</b>	FVC081822A																						
<b>Product Description:</b>	Lyophilized white solids containing 1 mg Select-HA (determined by Carbazole Assay)																						
<b>Quality Assay:</b>	<ul style="list-style-type: none"> <li>• Electrophoresis</li> <li>• Multi-angle Laser Light Scattering-Size Exclusion Chromatography (MALLS-SEC)</li> <li>• Chromogenic Limulus Amebocyte Lysate (LAL) Assay</li> </ul>																						
<b>Quality Parameters</b>	<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>1.01</td> <td>✓</td> </tr> <tr> <td>C. Molecular Mass</td> <td>360-440 kDa</td> <td>400 kDa</td> <td>✓</td> </tr> <tr> <td>D. Endotoxin Level</td> <td>≤0.1 EU/mg</td> <td>0.04 EU/mg</td> <td>✓</td> </tr> </tbody> </table>	Parameter	Required	Actual	Pass	A. Clarity on Gel	Clear	Clear	✓	B. Polydispersity	≤1.05	1.01	✓	C. Molecular Mass	360-440 kDa	400 kDa	✓	D. Endotoxin Level	≤0.1 EU/mg	0.04 EU/mg	✓	 <p><b>Electrophoresis Analysis</b> Select-HA was run on 1% (w/v) agarose gel and stained with 0.005% Stains-All in 50% ethanol.</p> <p><b>Lane 1</b> – Select-HA LoLadder (HYA-LOLAD-20, LL201812)</p> <p><b>Lane 2</b> – 400 kDa Select-HA (HYA-0400, lot# FVC081822A)</p>	
Parameter	Required	Actual	Pass																				
A. Clarity on Gel	Clear	Clear	✓																				
B. Polydispersity	≤1.05	1.01	✓																				
C. Molecular Mass	360-440 kDa	400 kDa	✓																				
D. Endotoxin Level	≤0.1 EU/mg	0.04 EU/mg	✓																				

Passes All Functional Specifications:



Date: 10/9/2023

Angie M. Branch, Biology Manager

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



Ph: 866-588-0455  
Fax: 801-588-0497

Echelon-inc.com