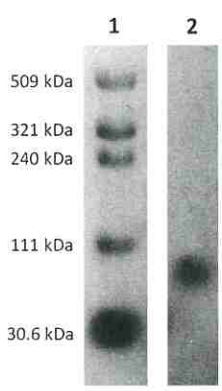


Echelon Biosciences Inc.

Certificate of Analysis

Product Name:	75 kDa Select-HA																						
Product Number:	HYA-0075																						
Lot Number:	FVC081822B																						
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>1.01</td> <td>✓</td> </tr> <tr> <td>C. Molecular Mass</td> <td>68-83 kDa</td> <td>74 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	68-83 kDa	74 kDa	✓	D. Endotoxin Level	≤0.1 EU/mg	0.04 EU/mg	✓	 <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 LoLadder (HYA-LOLAD-20, LL201812)</p> <p>Lane 2 – 75 kDa Select-HA (HYA-0075, lot# FVC081822B)</p>	
Parameter	Required	Actual	Pass																				
A. Clarity on Gel	Clear	Clear	✓																				
B. Polydispersity	≤1.05	1.01	✓																				
C. Molecular Mass	68-83 kDa	74 kDa	✓																				
D. Endotoxin Level	≤0.1 EU/mg	0.04 EU/mg	✓																				

Passes All Functional Specifications: ABC

Date: 2/8/2023

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