

Product Number:	103710
Lot Number:	019504
Expiration:	13-Jul-2022
Storage:	-80°C

**Laboratory Analysis**

Means of Analysis	Characteristic	Assumed* Concentration (µg/mL)	Measured Concentration (µg/mL)
HD-1 Analyzer	Antigen Concentration	1.00	0.756

\* Based on Vendor Certificate of analysis for raw antigen concentration.

Review/Approval

<u>Junp11 Park</u>	<u>QC Supervisor</u>	<u></u>	<u>18 Mar 2021</u>
Name	Title	Signature/Date	

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.