

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	1.00

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 12 Mar 2021
Signature/Date

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