

Product Number:	103710
Lot Number:	107306
Expiration:	14-Mar-2023
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Assumed* Concentration (ug/mL)	Measured Concentration (ug/mL)
HD-1 Analyzer	Antigen Concentration	1.00	0.868

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

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel

Signature/Date

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