

Product Number:	103373
Lot Number:	931112
Expiration:	7-Nov-2021
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Assumed* Concentration (mg/mL)	Measured Concentration (mg/mL)
HD-1 Analyzer	Antigen Concentration	1.00	1.27

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

Review/Approval

Shivani Goel

Name

QC Manager

Title

Shivani Goel 09 Oct 2020
Signature/Date

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