

Product Number:	103217
Lot Number:	026801
Expiration:	24-Sep-2022
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Measured Concentration (pg/mL)
HD-1 Analyzer	Calibrator Concentration	2184

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 27 Oct 2020
Signature/Date

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