

Product Number:	101734
Lot Number:	927002
Expiration:	27-Sep-2021
Storage:	-20°C

Laboratory Analysis

Means of Analysis	Characteristic	Measured Concentration (pg/mL)
HD-1 Analyzer	Calibrator Concentration	4991

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 26/01/2021
Signature/Date

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