


Product Number:	101734
Lot Number:	911207
Expiration:	22-Apr-2021
Storage:	-20°C

Laboratory Analysis

Means of Analysis	Characteristic	Measured Concentration (pg/mL)
HD-1 Analyzer	Calibrator Concentration	5847

Review/Approval

Junpil Park	QC Supervisor	 05 Feb 2021
Name	Title	Signature/Date

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