

Product Number:	102243
Lot Number:	006212
Expiration:	2-Mar-2022
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	218

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 31 Aug 2020

Name

Title

Signature/Date

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