

Product Number:	102243
Lot Number:	027611
Expiration:	2-Oct-2022
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	1419

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 03Jun2021

Signature/Date

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