

Product Number:	102260
Lot Number:	127303
Expiration:	30-Sep-2023
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	556

Review/Approval

Shivani Goel
Name

QC
Title

Shivani Goel 04Oct2021
Signature/Date

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