

<b>Product Number:</b>	102260
<b>Lot Number:</b>	105801
<b>Expiration:</b>	27-Feb-2023
<b>Storage:</b>	-80°C

**Laboratory Analysis**

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	693

Review/Approval

Shivani Goel  
~~Junpil Park~~

Manager  
Supervisor, QC

Shivani Goel

Name

Title

Signature/Date

SG 05 Aug 2021

SG 05 Aug 2021

05 Aug 2021

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