

<b>Product Number:</b>	100877
<b>Lot Number:</b>	016013
<b>Expiration:</b>	8-Jun-2022
<b>Storage:</b>	-80°C

**Laboratory Analysis**

Means of Analysis	Characteristic	Result (ug/mL)
HD-1 Analyzer	Concentration	49.9

Review/Approval

Junpil Park

Supervisor, QC



23 Feb 21

Name

Title

Signature/Date

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