

Product Number:	103216
Lot Number:	006211
Expiration:	2-Mar-2022
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-1 Analyzer	Concentration	426

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 31 Aug 2020.
Signature/Date

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