

Product Number:	102035
Lot Number:	119204
Expiration:	11-Jul-2023
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-1 Analyzer	Concentration	4.07

Review/Approval

Wenkai Tsai

Name

QC

Title



23 Jul 2021

Signature/Date

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