

Product Number:	103287
Lot Number:	506416
Expiration:	5-Mar-2031
Storage:	-80°C

Means of Analysis	Characteristic	Result (ug/mL)
HD-X Analyzer	Concentration	116

Review/Approval

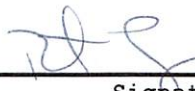
Robert Somogyi

QC Manager

Name

Title

Signature/Date

 26 MAR 2015

Quanterix Corporation

900 Middlesex Turnpike, Building 1, Billerica, MA 01821

techsupport@quanterix.com

©Quanterix Corp. Simoa® is a registered trademark of Quanterix Corp.

For research use only. Not for use in diagnostic procedures