

Product Number:	102260
Lot Number:	933806
Expiration:	4-Dec-2021
Storage:	-80°C

Laboratory Analysis

Means of Analysis	Characteristic	Result (ug/mL)
HD-1 Analyzer	Concentration	645

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler 24 Jul 2016

Name

Title

Signature/Date

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