

Product Number:	103705
Lot Number:	027201
Expiration:	28-Sep-2022
Storage:	-20°C

Laboratory Analysis

Concentration (pg/mL)	1223
------------------------------	------

Review/Approval

Shivani Goel

QC, Manager

Shivani Goel 03 Feb 2022

Name

Title

Signature/Date

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