

Product Number:	101734
Lot Number:	107906
Expiration:	20-Mar-2023
Storage:	-20°C

**Laboratory Analysis**

Concentration (pg/mL)	4973
-----------------------	------

Review/Approval

Shivani Goel

QC Manager

*Shivani Goel* 17 May 2022

Name

Title

Signature/Date

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