

Product Number:	102619
Lot Number:	026716
Expiration:	23-Sep-2022
Storage:	- 20°C

**Laboratory Analysis**

Concentration (pg/ml)	9343
-----------------------	------

Review/Approval Shivani Goel QC Shivani Goel 24 Aug 2021  
Name Title Signature/Date

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