

Product Number:	103217
Lot Number:	202410
Expiration:	24-Jan-2024
Storage:	-80°C

**Laboratory Analysis**

Concentration (pg/mL)	1417
-----------------------	------

Review/Approval

*Matt Smart*  
Name

QC  
Title

*MQA 3 May 2022*  
Signature/Date