

Product Number:	100895
Lot Number:	122105
Expiration:	9-Aug-2023
Storage:	- 80°C

Laboratory Analysis

Concentration (pg/ml)	3515
-----------------------	------

Review/Approval

Matt Smart Director, QC
Name Title

MAT 28Jun2022
Signature/Date