

Product Number:	101734
Lot Number:	130203
Expiration:	29-Oct-2023
Storage:	-20°C

Laboratory Analysis

Concentration (pg/mL)	4295
-----------------------	------

Review/Approval

Matt Koulouris

Name

QC Supervisor

Title

[Signature] 29 Mar 2023

Signature/Date