

Product Number:	102619
Lot Number:	220014
Expiration:	19-Jul-2024
Storage:	- 20°C

Laboratory Analysis

Concentration (pg/ml)	3887
-----------------------	------

Review/Approval

Matt Koulouris

Name

QC Supervisor

Title

[Signature] 29 Mar 2023

Signature/Date