

Product Number:	104102
Lot Number:	226601
Expiration:	23-Sep-2024
Storage:	-20°C

Laboratory Analysis

Concentration (pg/mL)	47,678
-----------------------	--------

Review/Approval

Tony Icm
Name

QL
Title


Signature/Date

14 FEB 2023