

Product Number:	100758 - 100765
Lot Number:	305311
Expiration:	22-Feb-2025
Storage:	-20°C

Calibrator Levels (pg/mL)	
A (100758)	0.00
B (100759)	0.077
C (100760)	0.246
D (100761)	0.873
E (100762)	2.61
F (100763)	8.43
G (100764)	25.8
H (100765)	64.4

Review/Approval Sonia Khillan QC Manager Sonia Khillan 14SEP2023  
Name Title Signature/Date