

Product Number:	103297
Lot Number:	231413
Expiration:	10-Nov-2023
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sonia Khillan QC Manager Sonia Khillan 11/06/2022  
Name Title Signature/Date