

Product Number:	103297
Lot Number:	303709
Expiration:	6-Feb-2024
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sania Khillan QC Manager Sania Khillan 08 FEB 2023
Name Title Signature/Date