

Product Number:	103297
Lot Number:	312405
Control Number:	312405A
Expiration:	4-May-2024
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 12 MAY 2023
 Name Title Signature/Date