

Product Number:	103297
Lot Number:	223506
Expiration:	23-Aug-2023
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sonia Khillan QC Sonia Khillan / 24 AUG 2022  
Name Title Signature/Date