

Product Number:	103297
Lot Number:	333109
Expiration:	27-Nov-2024
Storage:	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval Sonia Khillan Senior QC Analyst Sonia K.P./c 30 Nov 2023

Name Title Signature/Date