

<b>Product Number:</b>	103297
<b>Lot Number:</b>	108006
<b>Expiration:</b>	21-Mar-2022
<b>Storage:</b>	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

Shivani Goel 26 Mar 2021  
Signature/Date