

<b>Product Number:</b>	103297
<b>Lot Number:</b>	105514
<b>Expiration:</b>	24-Feb-2021
<b>Storage:</b>	2-8°C

Characteristic	
SBG Stock Concentration (nM)	10nM

Means of Analysis	Result
HD-1/ HDX	PASS

Review/Approval

Junpil Park

QC, Supervisor



11 Mar 2021

Name

Title

Signature/Date