

<b>Product Number:</b>	103397
<b>Lot Number:</b>	204911
<b>Expiration:</b>	8-Feb-2023
<b>Storage:</b>	2-8°C

**Laboratory Analysis**

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	50
---------------------------------------	----

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 28 Mar 2022  
Signature/Date