

Product Number:	103397
Lot Number:	124216
Expiration:	30-Aug-2022
Storage:	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* 18 May 2022  
Signature/Date