

Product Number:	103397
Lot Number:	124215
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  
Title

*Shivani Goel 30 Dec 2021*  
Signature/Date