

Product Number:	103397
Lot Number:	230622
Expiration:	29-Aug-2023
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

*Matt Koulouris*  
Name

*QC Supervisor*  
Title

*M. Koulouris* 07Feb2023  
Signature/Date