

Product Number:	103397
Lot Number:	307305
Expiration:	14-Feb-2024
Storage:	2-8°C

**Laboratory Analysis**

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval Brian Ciccariello VP QA+RA B-C 19 APR 2023  
Name Title Signature/Date