

Product Number:	103397
Lot Number:	222904
Expiration:	17-Aug-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval	<u> Matt Kalouris </u>	<u> QC supervisor </u>	<u> MK 02 Dec 2022 </u>
	Name	Title	Signature/Date