

Product Number:	104433
Lot Number:	315715
Control Number:	315715A
Expiration:	6-Jun-2026
Storage:	-20°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD-X Analyzer	PASS

SBG Concentrate Concentration (nM)	65
------------------------------------	----

Review/Approval

*Sonia Khillan*  
Name

*QC Manager*  
Title

*Sonia Khillan 10 OCT 2023*  
Signature/Date