

Product Number:	104433
Lot Number:	315714
Control Number:	315714A
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





Name

Title

Signature/Date