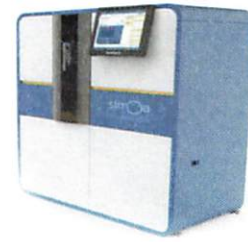


Product Number:	101735
Lot Number:	503872
Expiration:	14-Feb-2024



Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	310014
SBG Concentrate	103397	307305

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval

Ratish Suresh

Name

QC
Supervisor

Title

Ratish Suresh

Signature/Date

20 APR 2023