

Product Number:	101735
Lot Number:	503463
Expiration:	8-Feb-2023



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

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	211404
SBG Concentrate	103397	204912

Laboratory Analysis

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

Review/Approval

Richa Shrestha

QC Supervisor

Name

Title

27Apr2022

Signature/Date