

Product Number:	103615
Lot Number:	503970
Paired Reagent Kit	503866
Expiration:	18-Jul-2024
Platform(s):	HD-1 / HD-X & SR-X



Component	Part Number	Lot Number
N2PB Sample Diluent	103516	323612
Calibrator Diluent	103614	321506
NF-light Calibrator Concentrate	102619	220014
GFAP Calibrator Concentrate	103622	219910

Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Calibrator Concentrate Value Assignment (pg/mL)	
NF-Light	5027
GFAP	103950

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan 19SEP2023

Signature/Date