

Product Number:	103671
Lot Number:	503380
Paired Reagent Kit	503248
Expiration:	24-Nov-2022
Platform(s):	HD-1 / HD-X & SR-X



Component	Part Number	Lot Number
N4PE Plasma Sample Diluent	103659	132817
N4PE CSF Sample Diluent	103727	132818
Calibrator Diluent	103661	203713
Abeta 40 Calibrator Concentrate	101734	107906
Abeta 42 Calibrator Concentrate	103217	211005
GFAP Calibrator Concentrate	103622	027001
NF-light Calibrator Concentrate	102619	026716

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

Calibrator Concentrate Value Assignment (pg/mL)	
Abeta 40	4229
Abeta 42	2125
GFAP	149367
NF-Light	9343

Review/Approval

Matt Smart

QC director

M. A. 01/15/2022

Name

Title

Signature/Date