

Product Number:	103671
Lot Number:	502782
Paired Reagent Kit	502721
Expiration:	27-Sep-2021
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
N4PE Plasma Sample Diluent	103659	103312
N4PE CSF Sample Diluent	103727	103314
Calibrator Diluent	103661	103413
Abeta 40 Calibrator Concentrate	101734	927002
Abeta 42 Calibrator Concentrate	103217	026801
GFAP Calibrator Concentrate	103622	027001
NF-light Calibrator Concentrate	102619	026716

Calibrator Concentrate Value Assignment (pg/mL)	
Abeta 40	3592
Abeta 42	2184
GFAP	149367
NF-Light	9343

Review/Approval

Junpil Park

Name

QC Supervisor

Title

 01 Mar 2021

Signature/Date