

Product Number:	103615
Lot Number:	503144
Paired Reagent Kit	503094
Expiration:	17-Aug-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
N2PB Sample Diluent	103516	123011
Calibrator Diluent	103614	122908
NF-light Calibrator Concentrate	102619	026716
GFAP Calibrator Concentrate	103622	027001

Calibrator Concentrate Value Assignment (pg/mL)	
NF-Light	6683
GFAP	133346

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 04 May
Signature/Date 2022