

Quanterix

900 Middlesex Turnpike
Building 1
Billerica, MA 01821

tel:617.301.9400
fax:617.301.9401

www.quoterix.com

Assay Kit:	Lilly Neuro 4-Plex E (Bulk)
Product Number:	103971
Lot Number:	503474
Expiration:	24Nov2022
Platform(s):	HD-X

Component	Part Number	Lot Number
Bead Reagent	103653	134108
Detector Reagent	103655	214406
SBG Reagent	103657	132812
Plasma Sample Diluent	103659	132817
Calibrators	103667	211607
Control 0	103846	211604
Control 1	103668	211603
Control 2	103669	211603
RGP Reagent¹	103159	N/A

¹ RGP is not Kit Lot Specific

Reference Calibrators

Abeta 40 (pg/mL)	Conc. (pg/mL)	AEB*	AEB %CV*
A	0.000	0.002	19%
B	1.11	0.004	14%
C	2.11	0.007	7%
D	4.41	0.016	7%
E	9.33	0.053	1%
F	20.3	0.171	5%
G	40.0	0.474	1%
H	69.7	0.966	4%
Abeta 42 (pg/mL)	Conc. (pg/mL)	AEB*	AEB %CV*
A	0.000	0.004	5%
B	0.363	0.009	1%
C	0.721	0.023	3%
D	1.47	0.066	5%
E	2.97	0.241	2%
F	6.06	0.850	0%
G	12.0	2.50	3%
H	20.5	6.49	7%
GFAP (pg/mL)	Conc. (pg/mL)	AEB*	AEB %CV*
A	0.000	0.003	32%
B	5.23	0.018	4%
C	13.6	0.051	7%
D	47.2	0.156	4%
E	133	0.460	1%
F	483	1.63	6%
G	1607	4.44	5%
H	4884	11.6	12%
NF-Light (pg/mL)	Conc. (pg/mL)	AEB*	AEB %CV*
A	0.000	0.010	11%
B	0.444	0.024	4%
C	1.48	0.056	3%
D	4.64	0.158	4%
E	13.0	0.449	0%
F	44.4	1.40	6%
G	134	3.77	2%
H	389	10.5	12%

*Data from kit release run

Control Concentrations from 5 Runs

Control 0	Run 1 Conc. (pg/mL)	Run 2 Conc. (pg/mL)	Run 3 Conc. (pg/mL)	Run 4 Conc. (pg/mL)	Run 5 Conc. (pg/mL)	Mean Conc. (pg/mL)	Conc. %CV
Abeta 40	8.64	8.14	8.54	9.76	8.35	8.68	7%
Abeta 42	2.56	2.67	2.66	3.06	2.54	2.70	8%
GFAP	95.1	100.9	97.7	94.9	99.6	97.7	3%
NF-Light	10.6	10.3	11.6	11.1	11.0	10.9	5%

Control 1	Run 1 Conc. (pg/mL)	Run 2 Conc. (pg/mL)	Run 3 Conc. (pg/mL)	Run 4 Conc. (pg/mL)	Run 5 Conc. (pg/mL)	Mean Conc. (pg/mL)	Conc. %CV
Abeta 40	17.5	16.8	17.5	17.6	17.2	17.3	2%
Abeta 42	5.31	5.42	5.16	5.43	5.30	5.32	2%
GFAP	185	197	188	202	194	193	4%
NF-Light	20.9	21.3	22.3	21.3	22.7	21.7	3%

Control 2	Run 1 Conc. (pg/mL)	Run 2 Conc. (pg/mL)	Run 3 Conc. (pg/mL)	Run 4 Conc. (pg/mL)	Run 5 Conc. (pg/mL)	Mean Conc. (pg/mL)	Conc. %CV
Abeta 40	132	132	132	133	134	132	1%
Abeta 42	29.7	30.3	30.0	30.9	29.7	30.1	2%
GFAP	3930	3938	3763	3667	3690	3798	3%
NF-Light	464	458	430	438	440	446	3%

Panel Results from Kit Release

Panel	Abeta 40 % Bias to Reference	Abeta 42 % Bias to Reference	GFAP % Bias to Reference	NF-Light % Bias to Reference
Panel 1	10%	1%	9%	8%
Panel 2	1%	8%	3%	3%
Panel 3	10%	2%	9%	2%
Panel 4	N/A	3%	4%	1%
Panel 5	13%	5%	7%	3%
Panel 6	3%	11%	7%	4%
Panel 7	16%	4%	4%	0%
Panel 8	N/A	2%	3%	1%
Panel 9	16%	12%	4%	N/A
Panel 10	14%	10%	5%	0%
Panel 11	9%	5%	0%	6%
Panel 12	N/A	8%	1%	1%
Panel 13	3%	4%	5%	N/A
Panel 14	11%	6%	9%	9%
Panel 15	N/A	8%	12%	3%
Panel 16	N/A	10%	2%	4%
Panel 17	8%	4%	2%	N/A
Panel 18	0%	2%	1%	6%
Panel 19	4%	1%	7%	0%
Panel 20	N/A	3%	13%	7%

Review/
Approval

Matt Smart Director QC

MMA 07/24/2022

Name

Title

Signature/Date