



CERTIFICATE OF ANALYSIS

Page 1 of 3

COA-04690

Product Description: A β Peptide Panel 1 (6E10) V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K15200E-1, K15200E-2, K15200E-4, K15200G-1, K15200G-2, K15200G-4

Kit Lot Number: K0040254

Expiration Date: 31 MAY 2016

Acceptance Data:

Metric	Specifications	Assay			
		A β 40	A β 38	A β 42	
% Recovery Concentration	Calibrators	80% - 120%	Pass	Pass	Pass
	Controls	75% - 125%	Pass	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass
% Recovery CV	Calibrators	<20%	Pass	Pass	Pass
	Controls	<20%	Pass	Pass	Pass
	ULOQ	<20%	Pass	Pass	Pass
	LLOQ	<20%	Pass	Pass	Pass
Non-specific binding		<1.0%	Pass	Pass	Pass

Concentration of Materials Tested:

Units = pg/mL	Assay		
	A β 40	A β 38	A β 42
40X Calibrator ^a	599 000	427 000	53 300
Neurodegeneration Control 1 (A00C0216)	8886	4442	1496
Neurodegeneration Control 2 (A00C0217)	2988	1097	489
Neurodegeneration Control 3 (A00C0218)	521	234	186
CSF Control 1 ^b	6913	4540	1687
CSF Control 2 ^b	3671	5183	568
CSF Control 3 ^b	1897	645	256
ULOQ	7000	8475	1271
LLOQ	25.0	60.0	3.13

^a Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentration, please refer to the Standard Curve Data on page 2.

^b CSF Controls are multi-analyte controls used to verify assay performance and precision in Human CSF test matrices. Please see Product Insert for experimental data.



CERTIFICATE OF ANALYSIS

Standard Curve Data from Kit QC:

A β 40		
Conc., (pg/mL)	Average Signal	%CV
14 975	661 110	5.3
3744	205 799	3.8
936	36 446	4.2
234	3830	5.1
58.5	474	5.8
14.6	139	4.3
3.66	92	5.8
0	81	7.3

A β 38		
Conc., (pg/mL)	Average Signal	%CV
10 675	796 351	5.0
2669	122 100	5.2
667	12 213	3.7
167	1334	4.8
41.7	282	5.6
10.4	143	6.6
2.61	113	7.1
0	106	8.4

A β 42		
Conc., (pg/mL)	Average Signal	%CV
1333	835 817	3.8
333	195 777	3.6
83.3	31 410	4.2
20.8	3967	5.1
5.21	735	5.2
1.30	201	5.1
0.325	99	6.3
0	70	8.0

A β Peptide Panel 1 (6E10) V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
A β Peptide Panel 1 Plate	Z0055037	2-8°C	30 NOV 2016
Anti-A β 6E10 Antibody	D0040447	2-8°C	31 OCT 2016
A β 1-38 Peptide	A0040085	≤ -70°C	31 MAY 2016
A β 1-40 Peptide	A0040100	≤ -70°C	31 OCT 2016
A β 1-42 Peptide	A0040101	≤ -70°C	31 OCT 2016
Diluent 35	M0350019	2-8°C	31 OCT 2016
Diluent 100	Y0270200	2-8°C	30 NOV 2016
A β 40 Blocker	R0110006	2-8°C	30 NOV 2016
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals	Not Kit Specific	Room Temperature	N/A
Neurodegeneration Control 1	Not Kit Specific; See Separate Control Pack CoA		
Neurodegeneration Control 2			
Neurodegeneration Control 3			




CERTIFICATE OF ANALYSIS

Page 3 of 3

COA-04690

Statement: The above kit lot was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dana Short	Quality Control		22 SEP 2014