CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human Total Tau Kit

Kit Catalog Numbers: K151LA-Series

Kit Lot Number: K0080661

Expiration Date: 30 Apr 2019

Acceptance Data/ Percent Recovery Specifications:

	Calibrator		Neurodegeneration Control 1		Neurodegeneration Control 2		Neurodegeneration Control 3		
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu Total Tau	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	cv	Value	Recovery	CV
1	hu Total Tau	30.0	80% - 120%	< 20%	8000	80% - 120%	< 20%

Stock Concentration of Material Tested:

Spot	Name	Item	Conc	Units
1	hu Total Tau	20X Calibrator	412000	pg/mL
		Neurodegeneration Control 1	1300	pg/mL
		Neurodegeneration Control 2	381	pg/mL
		Neurodegeneration Control 3	137	pg/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

hu Total Tau - Spot 1					
Conc. pg/mL	ECL Signal	%CV			
20600	411,000	2.6			
6870	150,000	2.5			
2290	45,300	2.1			
763	11,900	2.1			
254	2,760	3.9			
84.8	698	3.8			
28.3	222	4.1			
0.00	68	6.3			

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0055745	2-8°C	30 Apr 2019
Anti-Total Tau Antibody	D0080830	2-8°C	30 Apr 2019
Tau Calibrator	A0080120	≤ -70°C	31 Jul 2019
Diluent 35	M0350027	2-8°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus				
Plate Seals	Not Kit Specific			
Wash Buffer (20X), 100mL	Not Kit Specific			
Neurodegeneration Control 1	Not Kit Specific			
Neurodegeneration Control 2	Not Kit Specific			
Neurodegeneration Control 3	Not Kit Specific			

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	Dawn Stollar	12 OCT 2017