

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: K0080687
Expiration Date: 31 May 2019

Acceptance Data/ Percent Recovery Specifications:

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

Spot	Name	LLOQ			ULOQ		
		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

CERTIFICATE OF ANALYSIS

Page 2 of 2

Standard Curve Data:

hu Total Tau - Spot 1		
Conc. pg/mL	ECL Signal	%CV
20600	368,000	4.5
6870	152,000	4.3
2290	47,700	4.6
763	12,500	2.8
254	3,030	2.9
84.8	762	3.8
28.3	270	3.7
0.00	79	5.3

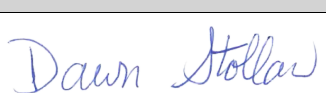
Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0055808	2-8°C	30 Jun 2019
Anti-Total Tau Antibody	D0080880	2-8°C	31 Jul 2019
Tau Calibrator	A0080120	≤ -70°C	31 Jul 2019
Diluent 35	M0350025	2-8°C	31 May 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		03 APR 2017