

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Human Total Tau Kit
Kit Catalog Numbers: K151LA-Series
Kit Lot Number: K0081086
Expiration Date: 31 Jul 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 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 Tau	30.0	80% - 120%	< 20%	8000	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Tau	20X Calibrator	412000	pg/mL
		Neurodegeneration Control 1	1340	pg/mL
		Neurodegeneration Control 2	406	pg/mL
		Neurodegeneration Control 3	171	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 Tau - Spot 1		
Conc. pg/mL	ECL Signal	%CV
20600	279,000	5.3
6870	108,000	5.1
2290	33,400	3.0
763	8,770	4.1
254	2,080	6.5
84.8	530	4.7
28.3	181	3.8
0.00	62	5.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0056362	2-8°C	31 Jan 2021
Anti-Total Tau Antibody	D0081077	2-8°C	31 Aug 2021
Tau Calibrator	A0080120	≤ -70°C	31 Jul 2019
Diluent 35	M0350029	2-8°C	30 Sep 2020
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	<i>Dawn Stollar</i>	11 SEP 2018