

**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:** K0081169

**Expiration Date:** 28 Feb 2021

**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	Tau	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	Tau	20X Calibrator	250000	pg/mL
		Neurodegeneration Control 1	1370	pg/mL
		Neurodegeneration Control 2	436	pg/mL
		Neurodegeneration Control 3	169	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:**

Tau - Spot 1		
Conc. pg/mL	ECL Signal	%CV
12500	176,000	6.4
4170	65,500	6.7
1390	19,000	4.9
463	4,770	6.6
154	1,100	6.0
51.4	296	6.7
17.1	123	4.9
0.00	69	9.1

## CERTIFICATE OF ANALYSIS

Page 2 of 2

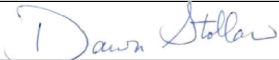
**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0056450	2-8°C	31 Jul 2021
Anti-Total Tau Antibody	D0081229	2-8°C	30 Jun 2022
Tau Calibrator	A0080179	≤ -70°C	28 Feb 2021
Diluent 35	M0350030	2-8°C	31 May 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100 mL	Not Kit Specific
Plate Seals	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
<b>Review/Approval</b>	Dawn Stollar	Quality		13 MAY 2019