

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: K0081302

Expiration Date: 31 Jan 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	189200	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
9460	178,000	3.8
3150	59,000	3.4
1050	15,600	3.0
350	3,560	4.4
117	884	4.9
38.9	291	4.2
13.0	140	5.4
0.00	80	15.0

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	D0081229	2-8°C	30 Jun 2022
Tau Calibrator	A0080196	≤ -70°C	31 Aug 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
Review/Approval	Karen Hamilla	Quality		19 NOV 2019