

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: K0081255

Expiration Date: 31 May 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	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	Total Tau	30.0	80% - 120%	< 20%	8000	80% - 120%	< 20%

Stock Concentration of Material Tested:

Spot	Name	Item	Conc	Units
1	Total 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:

Total Tau - Spot 1		
Conc. pg/mL	ECL Signal	%CV
9460	127,000	7.1
3150	41,600	5.3
1050	10,800	4.3
350	2,380	7.1
117	562	5.2
38.9	177	2.8
13.0	83	5.4
0.00	47	6.1

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0056520	2-8°C	31 Oct 2021
Anti-Total Tau Antibody	D0081309	2-8°C	31 Dec 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		25 JUL 2019