



# Meso Scale Diagnostics, L.L.C.

Description: V-PLEX® and V-PLEX Plus Human Total Tau Kit

Catalog Number(s): K151LAE/G-1/2/4

Lot #: K0040290

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of Lot # K0040290 has been extended to April 30, 2017.

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	18 FEB 2016

This memo supersedes all other statements of date of expiration for this kit, Lot # K0040290.

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

COA-06513
-----------

**Product Description:** V-PLEX® and V-PLEX Plus Human Total Tau Kit  
**Kit Catalog Numbers:** K151LAE-1, K151LAE-2, K151LAE-4, K151LAG-1, K151LAG-2, K151LAG-4  
**Kit Lot Number:** K0040290  
**Expiration Date:** 30 Apr 2017

**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	192000	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**

COA-06513
-----------

**Standard Curve Data:**

Tau - Spot 1		
Conc. pg/mL	ECL Signal	%CV
9600	229,000	9.6
3200	75,800	3.0
1070	19,200	3.4
356	4,140	4.7
119	817	5.4
39.5	202	5.2
13.2	84	5.9
0.00	46	15.0

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0055277	2-8°C	31 Jul 2017
Anti-Total Tau Antibody	D0040510	2-8°C	31 Aug 2017
Tau Calibrator	A0040097	≤ -70°C	30 Apr 2016
Diluent 35	M0350020	2-8°C	31 Dec 2016
MSD Read Buffer T (4X)	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 kit was tested using MSD procedures and it conforms to MSD's specifications. 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 Assurance	<i>Dawn Stollar</i>	18 FEB 2016