

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human Total Tau Kit
Kit Lot Number: K0081643
Expiration Date: 31 Mar 2022
Kit Catalog Numbers: K151LA-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Neurodegeneration Control 1 (4G8)		Neurodegeneration Control 2 (4G8)		Neurodegeneration Control 3 (4G8)	
		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 Calibrator:


Spot	Name	Conc.	Units
1	Tau	215000	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Total Tau Plate	Z0056864	2-8°C	30 Jun 2023
Anti-Total Tau Antibody	D0081538	2-8°C	29 Feb 2024
Tau Calibrator	A0080233	≤ -70°C	31 May 2022
Diluent 35	M0350031	2-8°C	31 Mar 2022

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		23 APR 2021