

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-22 Kit
Kit Catalog Numbers: K152WV-Series
Kit Lot Number: K0081050
Expiration Date: 31 Jul 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1(mouse) Control 1		TH17 Panel 1(mouse) Control 2		TH17 Panel 1(mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-22	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-22	1.58	75% - 125%	< 20%	380	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-22	Calibrator	573	pg/mL
		TH17 Panel 1(mouse) Control 1	423	pg/mL
		TH17 Panel 1(mouse) Control 2	134	pg/mL
		TH17 Panel 1(mouse) Control 3	41.7	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-22 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
573	311,000	3.3
143	67,000	2.7
35.8	15,100	2.4
8.95	3,720	1.5
2.24	1,060	2.9
0.560	366	1.9
0.140	196	2.0
0.00	136	2.2

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-22 Plate	Z0046871	2-8°C	31 Dec 2020
Anti-ms IL-22 Antibody	D0081049	2-8°C	31 Oct 2021
TH17 Panel 1(mouse) Calibrator Blend	A0080157	2-8°C	31 Oct 2022
Diluent 41	M0410022	≤ -10°C	31 Jul 2020
Diluent 45	M0450021	≤ -10°C	30 Sep 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
MSD Wash Buffer (20X), 100 mL	Not Kit Specific
TH17 Panel 1(mouse) Control 1	Not Kit Specific
TH17 Panel 1(mouse) Control 2	Not Kit Specific
TH17 Panel 1(mouse) 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	Dawn Stollar	Quality Assurance		09 AUG 2018