

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: K0080776

Expiration Date: 30 Sep 2019

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	555	pg/mL
		TH17 Panel 1 (mouse) Control 1	425	pg/mL
		TH17 Panel 1 (mouse) Control 2	136	pg/mL
		TH17 Panel 1 (mouse) Control 3	42.7	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
555	291,000	1.5
139	62,900	2.1
34.7	14,600	2.2
8.67	3,820	1.7
2.17	1,330	1.9
0.542	685	1.8
0.135	520	2.7
0.00	466	1.8

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-22 Plate	Z0046460	2-8°C	30 Sep 2019
Anti-ms IL-22 Antibody	D0080950	2-8°C	31 Oct 2019
TH17 Panel 1 (mouse) Calibrator Blend	A0080154	2-8°C	31 Oct 2022
Diluent 41	M0410018	≤ -10°C	30 Nov 2019
Diluent 45	M0450019	≤ -10°C	31 Dec 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
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	<i>Dawn Stollar</i>	02 JAN 2018