

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Liver Injury Panel 1 (rat) Kit
Kit Catalog Numbers: K15187D-series
Kit Lot Number: K0040406
Expiration Date: 31 Oct 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Liver Injury Panel 1 (rat) Control 1		Liver Injury Panel 1 (rat) Control 2		Liver Injury Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Arginase-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	alphaGST	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	Arginase-1	< 20.0%
4	alphaGST	< 1.0%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Arginase-1	20X Calibrator	4800	ng/mL
		Liver Injury Panel 1 (rat) Control 1	338	ng/mL
		Liver Injury Panel 1 (rat) Control 2	49.3	ng/mL
		Liver Injury Panel 1 (rat) Control 3	6.2	ng/mL
4	alphaGST	20X Calibrator	60200	ng/mL
		Liver Injury Panel 1 (rat) Control 1	4064	ng/mL
		Liver Injury Panel 1 (rat) Control 2	600	ng/mL
		Liver Injury Panel 1 (rat) Control 3	111	ng/mL

Control data was generated utilizing internal reference controls.

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

Page 2 of 2

Standard Curve Data:

Arginase-1 - Spot 1			alphaGST - Spot 4		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
240	510,000	3.1	3010	16,500	4.6
80.0	120,000	4.0	1000	13,600	4.8
26.7	28,000	4.5	334	9,160	4.3
8.89	7,920	7.0	111	4,760	5.1
2.96	2,530	6.5	37.2	2,190	5.0
0.988	932	4.3	12.4	994	6.6
0.329	404	4.4	4.13	547	6.1
0.00	161	6.5	0.00	296	6.5

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Liver Injury Panel 1 (rat) Plate	Z0056032	2-8°C	31 Jan 2020
Anti-rat Arginase-1 Antibody	D0040727	2-8°C	31 Jan 2020
Anti-rat alphaGST Antibody	D0040728	2-8°C	31 Mar 2020
Liver Injury Panel 1 (rat) Calibrator Blend	A0040201	≤ -70°C	31 Mar 2020
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
Diluent 100	Y0270249	2-8°C	31 Oct 2019
Blocker D-R	R0060022	≤ -10°C	31 Oct 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

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		07 MAR 2019