

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Liver Injury Panel 1 (rat) Kit

Kit Catalog Numbers: K15187D-Series

Kit Lot Number: K0040432

Expiration Date: 31 Mar 2020

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	Arg-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	alphaGST	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Arg-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:

Arg-1 - Spot 1			alphaGST - Spot 4		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
240	507,000	2.8	3010	14,200	2.5
80.0	142,000	4.3	1000	11,200	4.5
26.7	37,900	5.7	334	7,590	5.8
8.89	11,500	6.8	111	3,680	5.1
2.96	3,720	5.6	37.2	1,740	4.0
0.988	1,320	5.2	12.4	707	7.4
0.329	514	4.9	4.13	364	7.9
0.00	110	8.3	0.00	155	5.6

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Liver Injury Panel 1 (rat) Plate	Z0056291	2-8°C	30 Nov 2020
Anti-rat Arginase-1 Antibody	D0040778	2-8°C	30 Nov 2021
Anti-rat alphaGST Antibody	D0040779	2-8°C	30 Nov 2021
Liver Injury Panel 1 Calibrator Blend	A0040201	2-8°C	31 Mar 2020
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
Diluent 100	Y0270255	2-8°C	30 Sep 2020
Blocker D-R	R0010631	≤ -10°C	31 Oct 2020
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	<i>Dawn Stollar</i>	11 DEC 2018