

CERTIFICATE OF ANALYSIS

Product Description: Rat TIM-1/KIM-1/HAVCR Kit
Kit Catalog Numbers: K153JHC-Series
Kit Lot Number: K0040437
Expiration Date: 30 Nov 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Kidney Injury Panel (rat) Control 1		Kidney Injury Panel (rat) Control 2		Kidney Injury Panel (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Rat TIM-1/KIM-1/HAVCR	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Rat TIM-1/KIM-1/HAVCR	0.0200	80% - 120%	< 20%	10.0	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Rat TIM-1/KIM-1/HAVCR	20X Calibrator	192	ng/ml
		Kidney Injury Panel (rat) Control 1	19.9	ng/ml
		Kidney Injury Panel (rat) Control 2	3.3	ng/ml
		Kidney Injury Panel (rat) Control 3	0.972	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:

Rat TIM-1/KIM-1/HAVCR - Spot 1		
Conc. ng/ml	ECL Signal	%CV
9.59	682,000	1.9
3.20	249,000	4.6
1.07	75,400	3.7
0.355	23,700	3.9
0.118	7,860	4.6
0.0395	2,850	4.2
0.0132	1,220	2.5
0.00	451	2.5

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0056316	2-8°C	31 Dec 2020
Anti rat TIM-1/KIM-1/HAVCR Antibody	D0040792	2-8°C	31 Dec 2021
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040195	≤ -70°C	30 Nov 2019
Blocker A	Not Kit Specific	Not Kit Specific	N/A
Diluent 5	M0050064	≤ -10°C	28 Feb 2021
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		01 OCT 2018