

# CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** Rat TIM-1/KIM-1/HAVCR Kit

**Kit Catalog Numbers:** K153JHC-series

**Kit Lot Number:** K0040395

**Expiration Date:** 30 Apr 2019

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Kidney Injury Panel 1 (rat) Control 1		Kidney Injury Panel 1 (rat) Control 2		Kidney Injury Panel 1 (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 1 (rat) Control 1	20	ng/mL
		Kidney Injury Panel 1 (rat) Control 2	3.38	ng/mL
		Kidney Injury Panel 1 (rat) Control 3	0.712	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

**Standard Curve Data:**

Rat TIM-1/KIM-1/HAVCR - Spot 1		
Conc. ng/mL	ECL Signal	%CV
9.59	804,000	3.7
3.20	319,000	4.2
1.07	104,000	1.7
0.355	32,600	4.7
0.118	11,300	2.2
0.0395	3,970	2.5
0.0132	1,560	2.4
0.00	321	3.0


**CERTIFICATE OF ANALYSIS**

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0055951	2-8°C	31 Oct 2019
Anti-rat TIM-1/KIM-1/HAVCR Antibody	D0040703	2-8°C	30 Apr 2020
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040195	≤ -70°C	30 Nov 2019
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
Diluent 5	M0050060	≤ -10°C	30 Apr 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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		07 MAR 2019