

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** Rat TIM-1/KIM-1/HAVCR Kit  
**Kit Catalog Numbers:** K153JHC-Series  
**Kit Lot Number:** K0040427  
**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	173	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

**CERTIFICATE OF ANALYSIS**

**Standard Curve Data:**

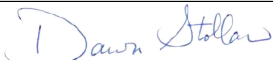
Rat TIM-1/KIM-1/HAVCR - Spot 1		
Conc. ng/mL	ECL Signal	%CV
8.64	935,000	1.8
2.88	349,000	1.9
0.960	109,000	2.8
0.320	34,200	2.3
0.107	11,600	1.9
0.0356	4,020	2.9
0.0119	1,540	2.4
0.00	310	3.7

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0056257	2-8°C	31 Oct 2020
Anti rat TIM-1/KIM-1/HAVCR Antibody	D0040766	2-8°C	30 Nov 2021
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040213	≤ -70°C	31 Oct 2020
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		06 MAR 2019