

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** Rat TIM-1/KIM-1/HAVCR Kit  
**Kit Catalog Numbers:** K153JHC-Series  
**Kit Lot Number:** K0040375  
**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

Controls tested may not be the same lot as shipped.

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
8.63	976,000	2.0
2.88	367,000	2.8
0.959	116,000	3.7
0.320	35,600	2.2
0.107	11,900	1.9
0.0355	4,160	2.4
0.0118	1,640	2.5
0.00	438	2.9

## CERTIFICATE OF ANALYSIS

Page 2 of 2

**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0055761	2-8°C	30 Apr 2019
Anti Rat TIM-1/KIM-1/HAVCR Antibody	D0040656	2-8°C	30 Apr 2019
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040186	≤ -70°C	31 May 2019
Blocker A Solution	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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 FEB 2017