

CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-07214

Product Description: Rat TIM-1/KIM-1/HAVCR Kit
Kit Catalog Numbers: K153JHC-1, K153JHC-2, K153JHC-4
Kit Lot Number: K0040333
Expiration Date: 31 Aug 2018

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Control 1		Control 2		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	213	ng/ml
		Control 1	20.0	ng/ml
		Control 2	3.38	ng/ml
		Control 3	0.710	ng/ml

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

COA-07214

Standard Curve Data:

Rat TIM-1/KIM-1/HAVCR - Spot 1		
Conc. ng/ml	ECL Signal	%CV
10.7	546,000	2.8
3.55	207,000	2.9
1.18	67,800	3.3
0.394	22,100	3.0
0.131	7,620	3.2
0.0438	2,780	3.6
0.0146	1,100	4.4
0.00	301	3.3

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0055586	2-8°C	31 Aug 2018
Anti-rat TIM-1/KIM-1/HAVCR Antibody	D0040596	2-8°C	31 Aug 2018
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040154	≤ -70°C	31 Aug 2018
Blocker A	Not Kit Specific	Room Temperature	N/A
Diluent 5	M0050058	≤ -10°C	30 Sep 2018
Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A

Statement:

The above kit was tested using MSD procedures and it conforms to MSD's specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	15AUG2016