



RFE MEMO

Page 1 of 1

Product Description(s): Rat TIM-1/KIM-1/HAVCR Kit

Catalog Number(s) / Size(s): K153JHC-Series

Lot Number: K0040466

Based on additional testing, the original date of expiration for this kit has been extended by 24 months. The expiration date of K0040466 has been extended to 30 APR 2023.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0056497	30 Sep 2021	30 APR 2023
Anti-rat TIM-1/KIM-1/HAVCR Antibody	D0040837	30 Sep 2022	30 APR 2023
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040223	30 Apr 2021	30 APR 2023
Diluent 5	M0050067	30 Nov 2021	30 APR 2023

Function	Name	Signature	Date
Quality	Karen Hamilla		02 MAY 2022

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Rat TIM-1/KIM-1/HAVCR Kit
Kit Catalog Numbers: K153JHC-Series
Kit Lot Number: K0040466
Expiration Date: 30 Apr 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TIM-1/KIM-1/HAVCR Control 1		TIM-1/KIM-1/HAVCR Control 2		TIM-1/KIM-1/HAVCR Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	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	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	TIM-1/KIM-1/HAVCR	20X Calibrator	180	ng/mL
		TIM-1/KIM-1/HAVCR Control 1	3	ng/mL
		TIM-1/KIM-1/HAVCR Control 2	0.45	ng/mL
		TIM-1/KIM-1/HAVCR Control 3	0.1	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:

TIM-1/KIM-1/HAVCR - Spot 1		
Conc. ng/mL	ECL Signal	%CV
9.00	843,000	2.3
3.00	319,000	2.8
1.00	100,000	3.1
0.333	31,200	3.5
0.111	10,300	3.9
0.0370	3,690	3.0
0.0123	1,470	3.8
0.00	408	2.9

CERTIFICATE OF ANALYSIS


Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0056497	2-8°C	30 Sep 2021
Anti-rat TIM-1/KIM-1/HAVCR Antibody	D0040837	2-8°C	30 Sep 2022
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040223	≤ -70°C	30 Apr 2021
Blocker A Kit	Not Kit Specific	Not Kit Specific	N/A
Diluent 5	M0050067	≤ -10°C	30 Nov 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	Karen Hamilla	Quality		21 NOV 2019