



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: K0040522

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0040522 has been extended to 31 JAN 2024.

The expiration date for the following kit component has been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Diluent 5	M0050070	31 Jan 2023	31 JAN 2024
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040255	28 Feb 2023	31 JAN 2024
Rat TIM-1/KIM-1/HAVCR Plate	Z0056769	30 Sep 2023	31 JAN 2024

Function	Name	Signature	Date
Quality	Karen Hamilla		17 NOV 2022

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

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Rat TIM-1/KIM-1/HAVCR Kit
Kit Lot Number: K0040522
Expiration Date: 31 Jan 2023
Kit Catalog Numbers: K153JHC-Series

Precision and Accuracy 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 Calibrator:


Spot	Name	Conc.	Units
1	TIM-1/KIM-1/HAVCR	192	ng/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat TIM-1/KIM-1/HAVCR Plate	Z0056769	2-8°C	30 Sep 2023
Rat TIM-1/KIM-1/HAVCR Antibody	D0040969	2-8°C	30 Nov 2025
Rat TIM-1/KIM-1/HAVCR Calibrator	A0040255	≤ -70°C	28 Feb 2023
Diluent 5	M0050070	≤ -10°C	31 Jan 2023

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		06 JUL 2022