



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Rat TIM-1/KIM-1/HAVCR Calibrator

**Catalog Number:** C03JH-2

**Lot Number:** A0040255

**Expiration Date:** 28 FEB 2023


**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
TIM-1/KIM-1/HAVCR	191.8	ng/mL

**Dilution to Top of Curve:** 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		30 OCT 2020