



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C03JH-2

Lot Number: A0040195

Expiration Date: 30 NOV 2019

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

Concentration of Materials Tested:

	Assay
	Rat TIM-1/KIM-1/HAVCR
Calibrator (ng/mL)	191.8

Statement: This 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>	18 OCT 2017