



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** Rat TIM-1/KIM-1/HAVCR Calibrator  
**Catalog Number:** C03JH-2  
**Lot Number:** A0040223  
**Expiration Date:** 30 APR 2021  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration of Materials Tested:**

	Assay
	<b>Rat TIM-1/KIM-1/HAVCR</b>
Calibrator ng/mL	180

Concentration: 20X

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

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 DEC 2018