



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Kidney Injury Panel 1 (rat) Calibrator Blend

**Catalog Number:** C0162-2

**Lot Number:** A0040225

**Expiration Date:** 30 APR 2021

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

**Concentration of Materials Tested:**

	Assay			
	Lipocalin-2	Osteopontin	Albumin	TIM-1
Calibrator (ng/mL)	4600	1036	980000	197.2

**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>	23 JAN 2019