



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** C0162-2

**Lot Number:** A0040240

**Expiration Date:** 30 APR 2022


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

### Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Lipocalin-2	4180	ng/mL
Osteopontin	928	ng/mL
Albumin	890000	ng/mL
TIM-1	179.4	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
<b>Review/Approval</b>	Karen Hamilla	Quality		28 JAN 2020