

**CERTIFICATE OF ANALYSIS**


**Product Description:** Kidney Injury Panel 3 (human) Calibrator Blend  
**Catalog Number:** C0189-2  
**Lot Number:** A0040295  
**Expiration Date:** 30 NOV 2027  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
Calbindin	538000	pg/mL
Clusterin	4160000	pg/mL
KIM-1	382000	pg/mL
Osteoactivin	758000	pg/mL
TFF3	35800	pg/mL
VEGF	43600	pg/mL

**Dilution to Top of Curve** 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Michelle Likiam	Quality		20 JUN 2025