

**CERTIFICATE OF ANALYSIS**

Product Description: Kidney Injury Panel 3 (human) Calibrator Blend
Catalog Number: C0189-2
Lot Number: A0040301
Expiration Date: 30 NOV 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Calbindin	536000	pg/mL
Clusterin	3840000	pg/mL
HAVCR1/KIM-1	370000	pg/mL
Osteoactivin	756000	pg/mL
TFF3	35200	pg/mL
VEGF-A	43200	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
Review/Approval	Karen Hamilla	Quality		05 AUG 2026