



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Kidney Injury Panel 3 (human) Calibrator Blend

Catalog Number: C0189-2

Lot Number: A0040271

Expiration Date: 31 DEC 2024

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Calbindin	504000	pg/mL
Clusterin	3780000	pg/mL
KIM-1	378000	pg/mL
Osteoactivin	736000	pg/mL
TFF3	37000	pg/mL
VEGF	46600	pg/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
Review/Approval	Karen Hamilla	Quality		22 MAR 2023