



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0189-2

Lot Number: A0040251

Expiration Date: 28 FEB 2023

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
α GST	662000	pg/ml
Calbindin	418000	pg/ml
Clusterin	3880000	pg/ml
KIM-1	450000	pg/ml
Osteoactivin	694000	pg/ml
TFF3	37200	pg/ml
VEGF	43600	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	Carissa B. Hubbard	Quality		13 NOV 2020