



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0189-2

Lot Number: A0040239

Expiration Date: 31 JUL 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
α GST	534000	pg/ml
Calbindin	428000	pg/ml
Clusterin	4000000	pg/ml
KIM-1	510000	pg/ml
Osteoactivin	736000	pg/ml
TFF3	48000	pg/ml
VEGF	47600	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		28 JAN 2020