



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0189-2

Lot Number: A0040212

Expiration Date: 31 JAN 2021

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
α GST	328000	pg/ml
Calbindin	644000	pg/ml
Clusterin	4960000	pg/ml
KIM-1	530000	pg/ml
Osteoactivin	956000	pg/ml
TFF3	48000	pg/ml
VEGF	56600	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	Xiomara Talavera	Quality		11 DEC 2019