



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Kidney Injury Panel 5 (human) Calibrator Blend

Catalog Number: C0188-2

Lot Number: A0040230

Expiration Date: 31 AUG 2022

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Albumin	4280000	pg/mL
B2M	434000	pg/mL
Cystatin C	738000	pg/mL
EGF	10380	pg/mL
NGAL	206000	pg/mL
OPN	2260000	pg/mL
UMOD	2300000	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		06 MAR 2020