



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: C0188-2

Lot Number: A0040210

Expiration Date: 31 AUG 2020

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

Concentration of Materials Tested:

	Assay						
	Albumin	B2M	Cystatin C	EGF	NGAL	OPN	UMOD
Calibrator (pg/mL)	3960000	416000	812000	10340	218000	1928000	2120000

Concentration: 20X

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 JUN 2018