



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07999

Product Description: Kidney Injury Panel 3 (human) Calibrator Blend
Catalog Number: C0189-2
Lot Number: A0040188
Expiration Date: 31 AUG 2019
Storage Temperature: ≤ -70°C

Concentration of Materials Tested:

	Assay						
	aGST	Calbindin	Clusterin	KIM-1	Osteoactivin	TFF3	VEGF
20X Calibrator (pg/mL)	38 000	498 000	4040 000	394 000	768 000	36 000	47 800

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 MAR 2017