

**CERTIFICATE OF ANALYSIS****Product Description:** Kidney Injury Panel 3 (human) Kit**Lot Number:** K0040581**Expiration Date:** 30 JUN 2027**Catalog Number:** K15189D-series**Lot-Specific Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Calbindin Antibody	D0041073	2-8°C	31 JUL 2028
Anti-hu Clusterin Antibody	D0041074	2-8°C	31 JUL 2028
Anti-hu Osteoactivin Antibody	D0041076	2-8°C	31 JUL 2028
Anti-hu TFF3 Antibody	D0041077	2-8°C	31 JUL 2028
Anti-hu VEGF Antibody	D0041078	2-8°C	31 JUL 2028
Human HAVCR1/KIM-1 Antibody	D0041075	2-8°C	31 AUG 2028
Kidney Injury Panel 3 (human) Plate	Z0057549	2-8°C	31 JUL 2027
Diluent 37	M0370041	≤ -10°C	30 JUN 2027
Kidney Injury Panel 3 (human) Calibrator Blend	A0040295	≤ -70°C	30 NOV 2027

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Kidney Injury Panel 3 Control 1		Kidney Injury Panel 3 Control 2	
		Recovery	CV	Recovery	CV	Recovery	CV
2	Calbindin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	KIM-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	Osteoactivin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	TFF3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
7	VEGF	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

Spot	Name	Spec
2	Calbindin	< 1.0 %
3	Clusterin	< 1.0 %
4	KIM-1	< 1.0 %
5	Osteoactivin	< 1.0 %
6	TFF3	< 1.0 %
7	VEGF	< 1.0 %

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
2	Calbindin	538000	pg/mL
3	Clusterin	4160000	pg/mL
4	KIM-1	382000	pg/mL
5	Osteoactivin	758000	pg/mL
6	TFF3	35800	pg/mL
7	VEGF	43600	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		16 JUL 2025