

**CERTIFICATE OF ANALYSIS**

Product Description: Kidney Injury Panel 3 (human) Kit
Lot Number: K0040572
Expiration Date: 31 MAR 2026
Catalog Number: K15189D-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Calbindin Antibody	D0041055	2-8°C	29 FEB 2028
Anti-hu Clusterin Antibody	D0041056	2-8°C	29 FEB 2028
Anti-hu Osteoactivin Antibody	D0041058	2-8°C	30 NOV 2027
Anti-hu TFF3 Antibody	D0041059	2-8°C	30 NOV 2027
Anti-hu VEGF Antibody	D0041060	2-8°C	30 NOV 2027
Human HAVCR1/KIM-1 Antibody	D0041057	2-8°C	30 NOV 2027
Kidney Injury Panel 3 (human) Plate	Z0057478	2-8°C	31 DEC 2026
Diluent 37	M0370040	≤ -10°C	31 OCT 2026
Kidney Injury Panel 3 (human) Calibrator Blend	A0040283	≤ -70°C	31 MAR 2026

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%

Non-Specific Binding Specifications:

Spot	Name	Spec
2	Calbindin	< 1.0%
3	Clusterin	< 1.0%
4	HAVCR1/KIM-1	< 1.0%
5	Osteoactivin	< 1.0%
6	TFF3	< 1.0%
7	VEGF-A	< 1.0%




CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
2	Calbindin	488000	pg/mL
3	Clusterin	3840000	pg/mL
4	KIM-1	382000	pg/mL
5	Osteoactivin	758000	pg/mL
6	TFF3	37800	pg/mL
7	VEGF	46400	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		01 OCT 2024