



CERTIFICATE OF ANALYSIS

Product Description: Kidney Injury Panel 3 (human) Kit
Lot Number: K0040588
Expiration Date: 30 JUN 2027
Catalog Number: K15756D-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu Calbindin Antibody	D0041085	2-8°C	31 MAR 2029
Anti-hu Clusterin Antibody	D0041086	2-8°C	28 FEB 2029
Anti-hu Osteoactivin Antibody	D0041088	2-8°C	28 FEB 2029
Anti-hu TFF3 Antibody	D0041089	2-8°C	28 FEB 2029
Anti-hu VEGF-A Antibody	D0041090	2-8°C	28 FEB 2029
Human HAVCR1/KIM-1 Antibody	D0041087	2-8°C	28 FEB 2029
Kidney Injury Panel 3 (human) 96-Well 10-Spot SECTOR Plate	Z0057609	2-8°C	29 FEB 2028
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:

		Calibrator		Kidney Injury Panel 3 Control 1		Kidney Injury Panel 3 Control 2	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV
1	Calbindin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Clusterin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	Osteoactivin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	HAVCR1/KIM-1	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
9	TFF3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
10	VEGF-A	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%



CERTIFICATE OF ANALYSIS

Non-Specific Binding Specifications:

Spot	Name	Spec
1	Calbindin	< 1.0%
2	Clusterin	< 1.0%
3	Osteoactivin	< 1.0%
8	HAVCR1/KIM-1	< 1.0%
9	TFF3	< 1.0%
10	VEGF-A	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Calbindin	538000	pg/mL
2	Clusterin	4160000	pg/mL
3	Osteoactivin	758000	pg/mL
8	HAVCR1/KIM-1	382000	pg/mL
9	TFF3	35800	pg/mL
10	VEGF-A	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		12 NOV 2025