

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Kidney Injury Panel 3 (human) Kit
Kit Lot Number: K0040546
Expiration Date: 30 Nov 2024
Kit Catalog Numbers: K15189D-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		KIP-3 Control 2		KIP3 In-process Control 1 (Cal-2)		KIP3 In-process Control 1 (Cal-2)	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
2	Calbindin	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%
3	Clusterin	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%
4	KIM-1	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%
5	Osteoactivin	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%
6	TFF3	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%
7	VEGF	80% - 120%	< 20%	75% - 125%	75% - 125%	75% - 125%	< 20%	75% - 125%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB 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	504000	pg/mL
3	Clusterin	3780000	pg/mL
4	KIM-1	378000	pg/mL
5	Osteoactivin	736000	pg/mL
6	TFF3	37000	pg/mL
7	VEGF	46600	pg/mL


CERTIFICATE OF ANALYSIS

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Kidney Injury Panel 3 (human) Plate	Z0057205	2-8°C	30 Nov 2024
Anti-hu Calbindin Antibody	D0040989	2-8°C	30 Nov 2025
Anti-hu Clusterin Antibody	D0040990	2-8°C	30 Nov 2025
Anti-hu KIM-1 Antibody	D0040991	2-8°C	30 Nov 2025
Anti-hu Osteoactivin Antibody	D0040992	2-8°C	30 Nov 2025
Anti-hu TFF3 Antibody	D0040993	2-8°C	30 Nov 2025
Anti-hu VEGF Antibody	D0040994	2-8°C	30 Nov 2025
Kidney Injury Panel 3 (human) Calibrator Blend	A0040271	≤ -70°C	31 Dec 2024
Diluent 37	M0370036	≤ -10°C	31 Jan 2025

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

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		07 FEB 2023