

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Kidney Injury Panel 1 (rat) Kit

Kit Lot Number: K0040565

Expiration Date: 31 Aug 2026

Kit Catalog Numbers: K15162C-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Kidney Injury Panel 1 (rat) Plate	Z0057441	2-8°C	31 AUG 2026
Anti-rat Lipocalin-2 Antibody	D0041039	2-8°C	31 AUG 2027
Anti-rat Osteopontin Antibody	D0041040	2-8°C	31 AUG 2027
Albumin Tracer	D0041038	2-8°C	31 AUG 2027
Rat HAVCR1/KIM-1 Antibody	D0041041	2-8°C	31 AUG 2027
Kidney Injury Panel 1 (rat) Calibrator Blend	A0040284	≤ -70°C	31 JUL 2026
Diluent 29	M0290035	≤ -10°C	31 AUG 2026

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

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Kidney Injury Panel 1 Control 1		Kidney Injury Panel 1 Control 2		Kidney Injury Panel 1 Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	Lipocalin-2	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	Osteopontin	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	Albumin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	TIM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	Lipocalin-2	3.67	70% - 130%	< 20%	150	80% - 120%	< 20%
2	Osteopontin	0.390	80% - 120%	< 20%	37.5	80% - 120%	< 20%
3	Albumin	781	80% - 120%	< 20%	33200	80% - 120%	< 20%
4	TIM-1	0.0200	80% - 120%	< 20%	1.88	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	Lipocalin-2	Osteopontin	Albumin	TIM-1
1	Lipocalin-2	NA	< 5.0%	< 1.0%	< 3.0%
2	Osteopontin	< 1.0%	NA	< 1.0%	< 1.0%
3	Albumin	< 1.0%	< 1.0%	NA	< 1.0%
4	TIM-1	< 1.0%	< 1.0%	< 1.0%	NA


CERTIFICATE OF ANALYSIS

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Lipocalin-2	3300	ng/mL
2	Osteopontin	1044	ng/mL
3	Albumin	984000	ng/mL
4	TIM-1	53.6	ng/mL

Dilution to Top of Curve = 20X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 APR 2024