

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: Kidney Injury Panel 1 (rat) Kit
Kit Lot Number: K0040501
Expiration Date: 30 Apr 2022
Kit Catalog Numbers: K15162C-Series

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	< 1.0%	< 1.0%	< 1.0%
2	Osteopontin	< 5.0%	NA	< 1.0%	< 1.0%
3	Albumin	< 1.0%	< 1.0%	NA	< 1.0%
4	TIM-1	< 1.0%	< 1.0%	< 1.0%	NA

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	Lipocalin-2	4180	ng/mL
2	Osteopontin	928	ng/mL
3	Albumin	890000	ng/mL
4	TIM-1	179	ng/mL

CERTIFICATE OF ANALYSIS


Page 2 of 2

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Kidney Injury Panel 1 (rat) Plate	Z0056762	2-8°C	31 Jan 2023
Anti-rat Lipocalin-2 Antibody	D0040926	2-8°C	31 Jan 2024
Anti-rat Osteopontin Antibody	D0040927	2-8°C	31 Jan 2024
Albumin Tracer	D0040925	2-8°C	30 Apr 2024
Anti-rat TIM-1/KIM-1/HAVCR Antibody	D0040928	2-8°C	31 Jan 2024
Kidney Injury Panel 1 (rat) Calibrator Blend	A0040240	≤ -70°C	30 Apr 2022
Diluent 29	M0290027	≤ -10°C	31 Dec 2022

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	Karen Hamilla	Quality		04 MAR 2021