



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Kidney Injury Panel 1 (rat) Calibrator Blend

Catalog Number: C0162-2

Lot Number: A0040269

Expiration Date: 31 OCT 2024

Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Lipocalin-2	3270	ng/mL
Osteopontin	897	ng/mL
Albumin	906000	ng/mL
TIM-1	51.6	ng/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		27 JUN 2022