



CERTIFICATE OF ANALYSIS

Product Description: Kidney Injury Panel 1 (rat) Calibrator Blend
Catalog Number: C0162-2
Lot Number: A0040284
Expiration Date: 31 JUL 2026
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Lipocalin-2	3300	ng/mL
Osteopontin	1044	ng/mL
Albumin	984000	ng/mL
TIM-1	53.6	ng/mL

Dilution to Top of Curve: 20 X

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		04 MAR 2024