



CERTIFICATE OF ANALYSIS

Product Description: Kidney Injury Panel 1 (rat) Calibrator Blend
Catalog Number: C0162-2
Lot Number: A0040292
Expiration Date: 30 JUN 2028
Storage Temperature: ≤ -70°C

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
NGAL/LCN2	4280	ng/mL
Osteopontin	976	ng/mL
Albumin	1052000	ng/mL
HAVCR1/KIM-1	53.8	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		24 FEB 2026