



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Inflammation Panel 1 (rat) Calibrator Blend

Catalog Number: C0179-2

Lot Number: A0040273

Expiration Date: 28 FEB 2025

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

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
Lipocalin-2	994	ng/mL
TSP-1	41600	ng/mL
TIMP-1	180.2	ng/mL
MCP-1	260	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		25 OCT 2022