



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Inflammation Panel 1 (rat) Calibrator Blend

Catalog Number: C0179-2

Lot Number: A0040220

Expiration Date: 31 MAR 2021

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

Concentration of Materials Tested:

| | Assay | | | |
|-----------------------|-------------|--------|--------|-------|
| | Lipocalin-2 | TSP-1 | TIMP-1 | MCP-1 |
| Calibrator (ng/mL) | 1072 | 43 600 | 186.4 | 266 |

Concentration: 20X

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 19 NOV 2018 |