



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Proinflammatory Panel 1 (human) Calibrator Blend

Catalog Number: C0049-2

Lot Number: A0080172

Expiration Date: 30 JUN 2023

Storage Temperature: 2-8°C

Concentration of Materials Tested:

| | Assay | | | | | | | | | |
|--|---------------------|--------------------|------------|------------|------------|------------|-------------|----------------|-------------|---------------------|
| | hu IFN- γ | hu IL-1 β | hu IL-2 | hu IL-4 | hu IL-6 | hu IL-8 | hu IL-10 | hu IL-12p70 | hu IL-13 | hu TNF- α |
| Reconstituted Calibrator (pg/mL) | 1430 | 561 | 1430 | 227 | 750 | 616 | 337 | 550 | 513 | 346 |

Reconstitution Volume: 1000 μ L

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> | 11 JUL 2018 |