



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06361

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080071
Expiration Date: 30 NOV 2020
Storage Temperature: 2-8°C

Concentration of Materials Tested:

| | Assay | | | | | | | | | |
|--|---------------------|--------------------|------------|------------|------------|------------|--------------|-------------|----------------|---------------------|
| | ms IFN- γ | ms IL-1 β | ms IL-2 | ms IL-4 | ms IL-5 | ms IL-6 | ms KC/GRO | ms IL-10 | ms IL-12p70 | ms TNF- α |
| Reconstituted Calibrator (pg/mL) | 1000 | 1740 | 2540 | 1710 | 987 | 5680 | 2110 | 3150 | 36 800 | 628 |

Reconstitution Volume: 1.0 mL

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. 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> | 18DEC2015 |