



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-04736

Product Description: Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0048-2
Lot Number: A0080043
Expiration Date: 31 JUL 2019
Storage Temperature: 2-8°C

Concentration of Materials Tested:

Table with 11 columns: Assay (ms IFN-gamma, ms IL-1beta, ms IL-2, ms IL-4, ms IL-5, ms IL-6, ms KC/GRO, ms IL-10, ms IL-12p70, ms TNF-alpha) and one row for Reconstituted Calibrator (pg/mL) with values: 890, 1500, 2180, 1500, 869, 4200, 1840, 2800, 34 600, 532

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.

Table with 5 columns: Name, Function, Signature, Date. Row 1: Review/Approval, Caitlin Hansen, Quality Control, [Signature], 09 SEP 2014