



CERTIFICATE OF ANALYSIS

COA-03658

Product Description: Proinflammatory Panel 1 (rat) Calibrator Blend
Catalog Number: C0044-2
Lot Number: A0080005
Expiration Date: 31 MAR 2018

Concentration of Materials Tested:

Table with 11 columns: Units = pg/mL, Assay (rt IFN-gamma, rt IL-1beta, rt IL-2, rt IL-4, rt IL-5, rt IL-6, rt KC/GRO, rt IL-10, rt IL-13, rt TNF-alpha) and one row for Reconstituted Calibrator with values: 4470, 11 400, 46 000, 854, 9610, 10 600, 2580, 20 900, 1340, 936

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 and one row for Review/Approval with values: Dana Short, Quality Control, [Signature], 10 APR 2014