



CERTIFICATE OF ANALYSIS

COA - 03509

**Product Description:** MSD<sup>®</sup> Human ProInflammatory I 4-Plex Calibrator Blend

**Catalog #/Size:** 3-3004-020205 / 1 Plate

**Lot Number:** A0031429

**Expiration Date:** 15 Months From Date of Receipt

**Statement:** This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Source (host/species): Recombinant protein derived from DNA sequence expressed in *E. coli*

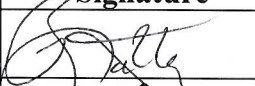

Concentration: 1 µg/mL

Purity: > 97% as determined by SDS-PAGE and visualized by silver stain.

Storage buffer: Diluent 1

Storage conditions: ≤ -70° C to -20° C

**The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.**

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		15 FEB 13
Review/Approval	Michael Eager	Quality Control		20 MAR 2013