



CERTIFICATE OF ANALYSIS

COA - 03511

Product Description: MSD[®] Human ProInflammatory 9-Plex Calibrator Blend

Catalog #/Size: C0007-2 / 1 Plate

Lot Number: A0031403

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): IL-1 β , IL-6, IL-8, TNF α , IFN γ , IL-2, GM-CSF: *E. coli*
IL-10, IL-12p70: *Sf21*

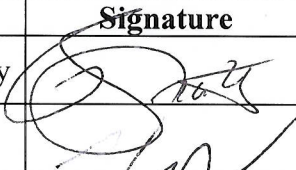

Concentration: 1 μ g/mL

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

Storage buffer: Diluent 1

Storage conditions: $\leq -70^{\circ}$ 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		20 FEB 13
Review/Approval	Michael Eager	Quality Control		05 APR 2013