

**CERTIFICATE OF ANALYSIS**

Product Description: Immune Response Panel 1 (human) Calibrator Blend
Catalog Number: C0742-2
Lot Number: A0080460
Expiration Date: 31 AUG 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
GM-CSF	6130	pg/mL
IFN- γ	8130	pg/mL
IL-1 β	3880	pg/mL
IL-2	14200	pg/mL
IL-6	5380	pg/mL
IL-12p70	9100	pg/mL
IL-10	5200	pg/mL
IL-8	4290	pg/mL
IL-17A	55500	pg/mL
TNF- α	12100	pg/mL

Dilution to Top of Curve: 5 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		19 JUN 2026