

**CERTIFICATE OF ANALYSIS**


Product Description: Immune Response Panel 1 (human) Calibrator Blend
Catalog Number: C0742-2
Lot Number: A0080422
Expiration Date: 30 SEP 2027
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
GM-CSF	7100	pg/mL
IFN- γ	7750	pg/mL
IL-1 β	3745	pg/mL
IL-2	14500	pg/mL
IL-6	5300	pg/mL
IL-12p70	9000	pg/mL
IL-10	5450	pg/mL
IL-8	4340	pg/mL
IL-17A	55000	pg/mL
TNF- α	12050	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	Michelle Likiam	Quality		16 JUN 2025