



CERTIFICATE OF ANALYSIS


Product Description: S-PLEX Proinflammatory Panel 1 (mouse) Calibrator Blend
Catalog Number: C0744-2
Lot Number: A00S0142
Expiration Date: 29 FEB 2028
Storage Temperature: $\leq -70^{\circ}\text{C}$

Concentration Assignment:

Analyte	Stock Concentration	Unit of Measure
IFN- γ	464000	fg/mL
IL-1 β	1496000	fg/mL
IL-2	5540000	fg/mL
IL-4	1072000	fg/mL
IL-5	3100000	fg/mL
IL-6	8500000	fg/mL
KC/GRO	1696000	fg/mL
IL-10	5540000	fg/mL
IL-12p70	68600000	fg/mL
TNF- α	2600000	fg/mL

Dilution to Top of Curve: 20 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 SEP 2025