



### CERTIFICATE OF ANALYSIS

**Product Description:** Cytolytic Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0741-2  
**Lot Number:** A0080420  
**Expiration Date:** 31 JUL 2027  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
GM-CSF	22200	pg/mL
Granzyme B	36400	pg/mL
IFN-γ	29700	pg/mL
IL-2	45700	pg/mL
IL-6	22700	pg/mL
Perforin	2280000	pg/mL
TNF-α	37200	pg/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
<b>Review/Approval</b>	Karen Hamilla	Quality		24 APR 2025