



**CERTIFICATE OF ANALYSIS**

**Product Description:** Cytolytic Panel 1 (human) Calibrator Blend  
**Catalog Number:** C0741-2  
**Lot Number:** A0080419  
**Expiration Date:** 31 JUL 2027  
**Storage Temperature:** ≤ -70°C

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
GM-CSF	22600	pg/mL
Granzyme B	37000	pg/mL
IFN- $\gamma$	29300	pg/mL
IL-2	47500	pg/mL
IL-6	22100	pg/mL
Perforin	2400000	pg/mL
TNF- $\alpha$	40100	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		25 MAR 2025