



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-06794

Product Description: Cytokine Panel 1 (human) Calibrator Blend
Catalog Number: C0050-2
Lot Number: A0080073
Expiration Date: 31 JAN 2021
Storage Temperature: 2-8°C

Concentration of Materials Tested:

	Assay									
	hu GM-CSF	hu IL-1 α	hu IL-5	hu IL-7	hu IL-12/23 p40	hu IL-15	hu IL-16	hu IL-17A	hu TNF- β	hu VEGF
Reconstituted Calibrator (pg/mL)	1130	487	883	972	3710	796	3400	6560	732	1260

Reconstitution Volume: 1.0 mL

Statement: The above product was manufactured and tested using MSD procedures and it conforms to MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	14 Apr 2016