



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Cytokine Panel 2 (human) Control Pack

Catalog Number: C4084-1

Lot Number: CP000052

Expiration Date: 31 JAN 2022

Concentration of Materials Tested:

Units = pg/mL	Assay							
	hu IL-17A/F	hu IL-17B	hu IL-17C	hu IL-17D	hu IL-1RA	hu IL-3	hu IL-9	hu TSLP
Reconstituted Control 1 (A00C0467)	5 913	487	2 360	5 479	670	3 173	1 272	581
Reconstituted Control 2 (A00C0468)	796	71	295	559	143	703	345	87
Reconstituted Control 3 (A00C0482)	74	133	77	146	50	240	66	12

Reconstitution Volume: 250 µL

Control concentrations were assigned across three kit lots.

Cytokine Panel 2 (human) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Control 1	A00C0467	2-8°C	31 JAN 2022
Cytokine Panel 2 (human) Control 2	A00C0468	2-8°C	31 JAN 2022
Cytokine Panel 2 (human) Control 3	A00C0482	2-8°C	28 FEB 2022

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>	25 JUL 2017