



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Proinflammatory Panel 1 (mouse) Control Pack

**Catalog Number:** C4048-1

**Lot Number:** CP000048

**Expiration Date:** 31 DEC 2021

### Concentration of Materials Tested:

Units = pg/mL	Assay									
	ms IFN- $\gamma$	ms IL-1 $\beta$	ms IL-2	ms IL-4	ms IL-5	ms IL-6	ms KC/GRO	ms IL-10	ms IL-12p70	ms TNF- $\alpha$
Reconstituted Control 1 (A00C0445)	282	663	951	437	322	1680	661	1030	13 800	152
Reconstituted Control 2 (A00C0446)	21.5	47.9	66.7	59.9	49.0	244	108	314	3930	29.1
Reconstituted Control 3 (A00C0447)	1.76	4.18	5.33	12.1	8.74	37.5	17.1	94.9	1080	5.10

**Reconstitution Volume:** 250  $\mu$ L

Proinflammatory Panel 1 (mouse) Control Pack Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (mouse) Control 1	A00C0445	2-8° C	31 DEC 2021
Proinflammatory Panel 1 (mouse) Control 2	A00C0446	2-8° C	31 DEC 2021
Proinflammatory Panel 1 (mouse) Control 3	A00C0447	2-8° C	31 DEC 2021

**Statement:** This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	07 MAR 2018