

BD Cytometric Bead Array

BD CBA Human and Mouse Th1/Th2/Th17 Kits

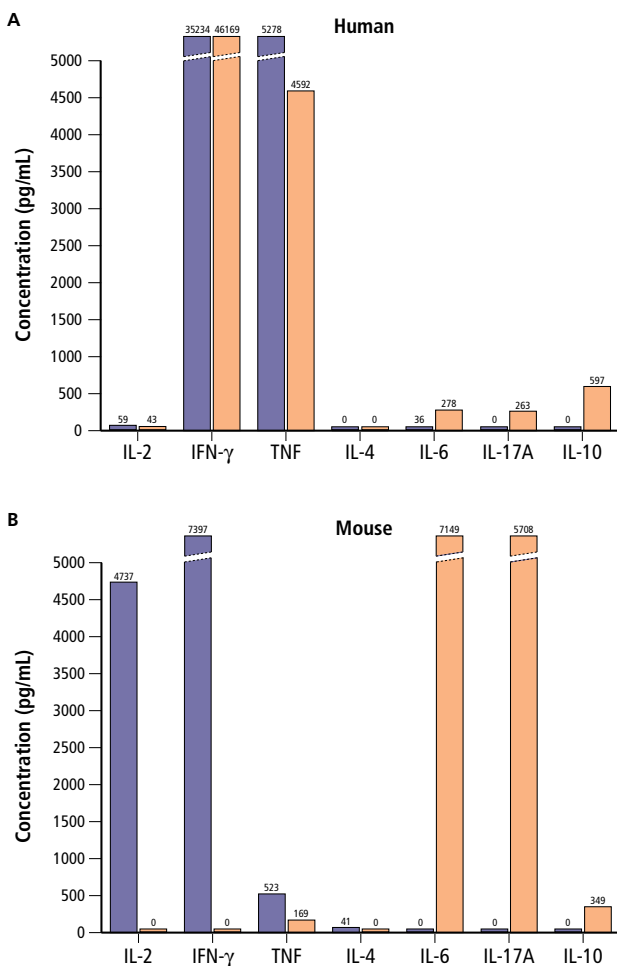
Features

Simultaneously quantitate IL-2, IL-4, IL-6, IL-10, IL-17A, IFN- γ , and TNF

Preconfigured for consistent results

Generate data from a single sample within hours

Compatible with most BD flow cytometers including those that have only a 488-nm laser



Measurement of Th17 cultures using the BD CBA Th1/Th2/Th17 Kits.

A) CD4⁺ human memory T cells isolated from whole blood were stimulated with plate-bound anti-CD3 and soluble anti-CD28 alone (blue) or in the presence of recombinant IL-23 (orange). **B)** Naïve CD4⁺ splenocytes were stimulated with plate-bound anti-CD3 and soluble anti-CD28 alone (blue) or in the presence of recombinant TGF- β and IL-6 (orange).

BD™ CBA Human and Mouse Th1/Th2/Th17 kits are essential tools for research in T cell biology. These kits contain reagents to measure particular sets of cytokines so that researchers can determine if T cell differentiation is being polarized toward a Th1, Th2, or Th17 response.

Cytokines measured

Each kit contains reagents to measure IL-2, IL-4, IL-6, IL-10, IL-17A, IFN- γ , and TNF. Many of these cytokines are multifunctional and are involved in proliferation (eg IL-2) and regulation (eg IL-10) of multiple T cell differentiation pathways. The kits feature key cytokines involved in differentiation of each cell type as well as the cytokines secreted by each cell type.

	Differentiation	Secretion
Th1	IFN- γ	IFN- γ , TNF
Th2	IL-4	IL-4, IL-6
Th17	IL-6	IL-17A, TNF

More data from less sample

The BD CBA system makes it possible to quantitate multiple analytes in serum, plasma, supernatant, and other complex biological matrices. The multiplex format significantly reduces sample requirements and experiment time. Multiplexing can also yield contextual answers that might be more relevant than single measurements.

The assay uses antibody-coated beads to capture analytes efficiently. Each bead has a unique fluorescence intensity that resolves as a unique population on a flow cytometer. Analyte bound to a bead is detected by a second antibody with a phycoerythrin (PE) label. The PE signal is proportional to the amount of bound analyte. Analysis software identifies each bead population, generates a standard curve, and calculates the concentration of unknown samples.

Ordering Information

Description	Cat. No.
Human Th1/Th2/Th17 CBA Kit	560484
Mouse Th1/Th2/Th17 CBA Kit	560485

Analysis Software

Description	Cat. No.
FCAP Array™ v1.0.2 (Mac OS®)	645447
FCAP Array™ v1.0.1 (PC Windows)	641488

Visit bdbiosciences.com/cba for more information.



BD Biosciences

2350 Qume Drive
San Jose, CA 95131
US Orders: 877.232.8995
bdbiosciences.com