

# Plasma Sample Diluent

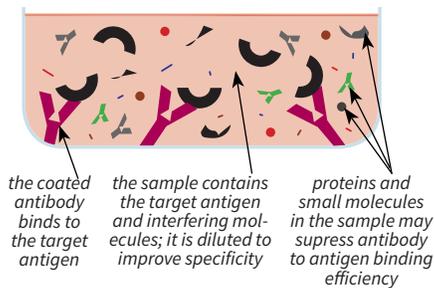
## A stable protein matrix for dilution of plasma and other aqueous-based biological samples.

Plasma Sample Diluent is formulated to provide a stable protein-friendly environment for dilution of neat plasma or serum samples for analysis in antibody-sandwich or antigen-down format ELISAs. It was specifically developed to provide a reliable solution to the occurrence of unwanted clotting events in plasma and serum samples within the microtiter plate wells. When assay sensitivity is a large concern, serum or plasma samples must be added to the ELISA plate neat or diluted only slightly, making clotting more likely. This issue can be essentially eliminated by diluting the potentially problematic samples 1:2 in Plasma Sample Diluent.

Sample diluents are used to dilute samples into the functional range of the assay and to create the standard curve. Due to the finite binding capacity of the plate-coated proteins, highly concentrated samples must be diluted in order to obtain absorbance readings within the sensitivity detection limits of the instrument. Properly formulated sample diluents will also reduce background noise associated with non-specific bridging of signal-generating conjugates to the microtiter plate surface.

Plasma Sample Diluent provides a BSA protein-buffered, neutral pH environment that is highly compatible with antibody-antigen interactions. Inclusion of a protein component helps minimize the degree of non-specific IgG adsorption onto coated and blocked ELISA plate wells during antigen-down screening of serum or plasma samples. Antimicrobial agents allow for room temperature bench-top use and extensive storage stability at 2-8°C.

### Dilute samples within the detection limits of the ELISA



### PLASMA SAMPLE DILUENT

Size	Catalog #
100 mL	#694
500 mL	#695
1 L	#696
10 L	#697

### INSTRUCTIONS:

- Generally, serum samples should be diluted at least 1:50 to minimize background noise caused by non-specific antibody binding.
- To dilute the sample 1:100, add 1 part sample to 99 parts Plasma Sample Diluent. For example, add 10  $\mu\text{L}$  of the sample to 990  $\mu\text{L}$  Plasma Sample Diluent for a total of 1,000  $\mu\text{L}$ .
- Highly concentrated samples may need to be diluted 1:1,000 or more.
- Once diluted, run the assay according to the specific ELISA protocol.
- Analyze the data. If the sample was diluted 1:100, the dilution factor must be considered when calculating the value. For example, if the sample generated an OD value that correlates to 500  $\text{pg/mL}$  based on the standard curve, multiply by the dilution factor of 100 to yield a true concentration of 50,000  $\text{pg/mL} = 50 \text{ ng/mL}$  in the sample.

For more ELISA protocols and information, please visit [www.immunochemistry.com](http://www.immunochemistry.com).

### SPECIFICATIONS:

- Clear to light yellow liquid
- 1X ready to use
- pH 7.2-7.6

### STORAGE:

- 24 months at 2-8°C
- 1 week at room temperature

### SAFETY & USAGE:

- Warning! May cause an allergic skin reaction.
- SDS available at [immunochemistry.com](http://immunochemistry.com)
- Not for human or drug use
- For research use only

*Build a better assay with ELISA Solutions from ImmunoChemistry Technologies.*



### BRIGHT MINDS, BRIGHT SOLUTIONS.

ImmunoChemistry Technologies, LLC gratefully acknowledges the significant contributions made by one of its founders, Brian W. Lee, Ph.D in the development of this product, including the creation and illustration of its strategy and protocol.

9401 James Avenue S., #155, Bloomington, MN 55431 USA

TOLL-FREE 1-800-829-3194 LOCAL 952-888-8788

FAX 952-888-8988

[WWW.IMMUNOCHEMISTRY.COM](http://WWW.IMMUNOCHEMISTRY.COM)