

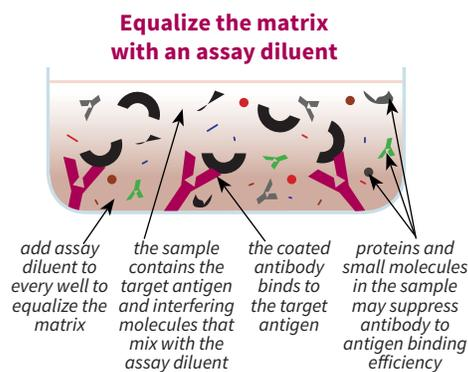
General Assay Diluent

Equalizes the sample and standard matrices for a more accurate result.

General Assay Diluent is formulated for testing serum, plasma, urine, and cell culture samples in all sandwich ELISA formats. Use of an assay diluent helps minimize matrix complexity differences between the sample (e.g., serum, etc.) and the diluent used to generate the standard curve of the ELISA.

Large differences between the sample and the standard diluent matrices will result in under-recovery of the target analyte present in sample wells. Complex and concentrated protein environments (the matrix) present in serum or plasma samples will greatly reduce the antigen-binding efficiency of the plate-adsorbed antibodies, resulting in a gross underestimation of the amount of target analyte present in the test samples. General Assay Diluent helps equalize the antibody-binding efficiencies between the standard curve and the sample wells.

To use, simply add 50-100 μ L to every well of the ELISA plate, including all wells designated for standards, controls, and samples. Then add the standards, controls, and samples to the plate. Mammalian protein additives included in the formulation serve to reduce non-specific interactions between the sample matrix proteins and the plate surface, thereby minimizing background noise. General Assay Diluent also inhibits complement and thrombin activity present in serum and plasma samples. Incorporation of an antimicrobial agent allows for room temperature bench-top use and extensive storage stability at 2-8°C. This assay diluent formulation ensures a more accurate and consistent ELISA performance, regardless of sample type.



GENERAL ASSAY DILUENT

Size	Catalog #
100 mL	#620
500 mL	#621
1 L	#622
10 L	#671

INSTRUCTIONS:

1. Dilute the standard curve, controls, and the samples as necessary. ICT offers several formulations of sample diluents in which to prepare the samples.
2. Pipette 50-100 μ L General Assay Diluent per well into every well of the plate.
3. Pipette 50-200 μ L of each standard, control, and sample into the plate.
4. Run the assay according to the specific ELISA protocol.
5. Analyze the data. Because all of the wells, including the standards and controls, received the same volume of assay diluent, there is no need to account for this dilution when calculating the results.

For more ELISA protocols and information, please visit 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
- 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