

Antigen-Down Assay Diluent

Minimizes clotting potential and reduces non-specific binding.

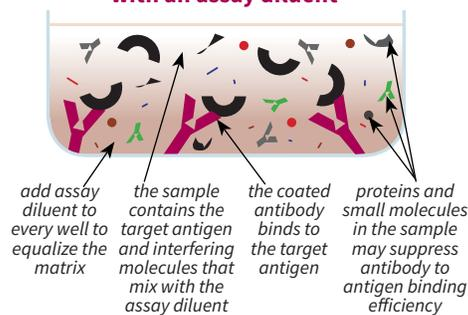
Antigen-Down Assay Diluent is formulated for testing serum and plasma samples in situations where an undesirable clotting event could occur while performing an antigen-down ELISA. Compared to ICT's other assay diluents, Antigen-Down Assay Diluent contains a higher molar concentration of calcium chelating agents and proprietary additives to eliminate the risk of clotting events in plasma samples and to minimize interference from complement and thrombin activity.

Antigen-Down Assay Diluent also provides a consistent protein matrix environment that will stabilize the antigen-specific target IgG analyte during high dilution ELISA titration protocols. Assessment of high titer, humoral IgG responses often requires dilutions >100,000-fold. This extensive dilution is often necessary to obtain a chromogen absorbance value that is within the sensitivity detection limits of the ELISA plate reader.

This novel assay diluent formulation includes mammalian protein and detergent additives to reduce the non-specific adsorption of generic serum IgG molecules onto the coated and blocked ELISA well surface. This feature can be especially critical when assessing sera derived from human, porcine, or bovine sources, which show a higher tendency to non-specifically adsorb. By minimizing such non-specific binding events, Antigen-Down Assay Diluent inhibits undesirable background noise and increases the sensitivity of the assay.

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. Incorporation of an antimicrobial agent into the formulation allows for room temperature bench-top use and extensive storage stability at 2-8°C.

Equalize the matrix with an assay diluent



ANTIGEN-DOWN ASSAY DILUENT

Size	Catalog#
100 mL	629
500 mL	630
1 L	631
10 L	674

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 Antigen-Down 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