

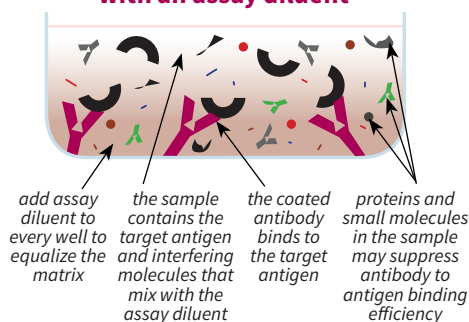
Neptune Assay Diluent

Provides a complex matrix for equalization of sample and standard matrices.

Neptune Assay Diluent is a non-mammalian protein-based reagent formulated to equalize complex matrix differentials that are often encountered when evaluating serum and plasma samples in antibody-sandwich ELISA formats. 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. Neptune Assay Diluent minimizes this serious issue in ELISA-format assays testing serum or plasma samples.

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. The diverse protein composition of Neptune Assay Diluent will minimize non-specific binding issues, decreasing background signal and increasing assay sensitivity. Neptune 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 novel assay diluent was designed to ensure more accurate and consistent ELISA performance, regardless of sample type.

Equalize the matrix with an assay diluent



NEPTUNE ASSAY DILUENT

Size	Catalog #
100 mL	#626
500 mL	#627
1 L	#628
10 L	#673

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 Neptune 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 rest of the assay according to the specific ELISA protocol.
5. Analyze the data. Because all 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:

- 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:

- Contains \leq 0.1% sodium azide
- SDS available at immunochemistry.com
- Product intended for research use or for further manufacturing into in vitro diagnostics reagents only.
- Not intended for use in human or therapeutics purposes.

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