

# Neptune™ HRP Conjugate Stabilizer, Non-Mammalian, 1X

Preserves the activity of HRP conjugates and reduces background signal.

Neptune HRP Conjugate Stabilizer is used to preserve concentrated stock conjugates, reconstitute lyophilized HRP conjugates, and dilute protein-peroxidase conjugates to their useful working titer in ELISAs and other immunology-based techniques.

Neptune HRP Conjugate Stabilizer, Non-Mammalian, 1X preserves the functional integrity of both the horseradish peroxidase enzyme and IgG components of the HRP-IgG conjugate complex. This proprietary formulation helps preserve the native three-dimensional conjugate structure during storage while concurrently enhancing the steric interaction or binding relationship between the IgG component of the HRP-IgG conjugate and the target antigen or antibody during the assay. By maintaining conjugate component activity and preventing contamination issues through the inclusion of an antimicrobial agent, Neptune HRP Conjugate Stabilizer, Non-Mammalian, 1X extends the functional utility of stored and reconstituted peroxidase conjugates. Extended stability studies clearly demonstrate this stabilizer's ability to preserve HRP-labeled IgG performance during long-term storage (see page 2).

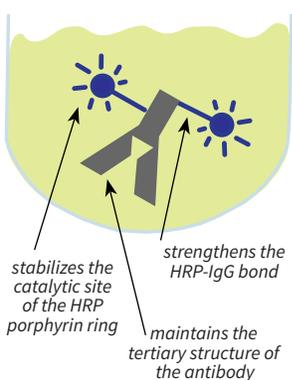
Neptune HRP Conjugate Stabilizer, Non-Mammalian, 1X is especially suited for ELISAs using anti-IgG HRP conjugates within traditional antigen-down or antibody sandwich formats. Incorporation of hydrolyzed non-mammalian proteins in place of traditional BSA protein

*continued on page 2...*

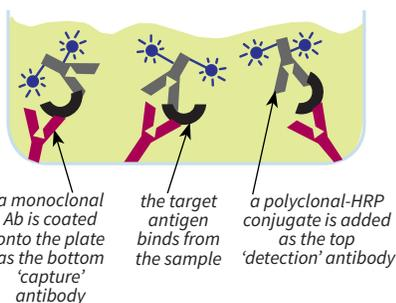
## BRIGHT MINDS, BRIGHT SOLUTIONS.

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.

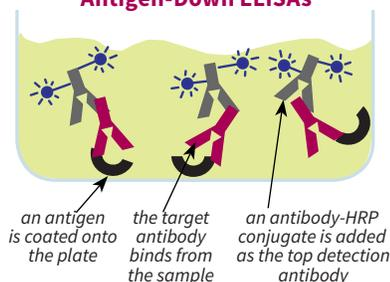
### Stabilizes HRP Conjugates



### Minimizes Background in Mono-Poly Sandwich ELISAs



### Minimizes Background in Antigen-Down ELISAs



## NEPTUNE™ HRP CONJUGATE STABILIZER, NON-MAMMALIAN, 1X

Size	Catalog #
100 mL	#6347
1 L	#6348
10 L	#6349

### INSTRUCTIONS:

1. Protect the stabilizer from light.
2. Gently mix the stabilizer; avoid bubbles.
3. Prepare the HRP conjugate solution at the predetermined dilution factor or concentration. For example, if preparing 50 mL at a dilution of 1:5,000, add 10 µL HRP-IgG conjugate stock concentrate to 49.99 mL Neptune HRP Conjugate Stabilizer, Non-Mammalian, 1X.
4. Mix for 15 minutes.
5. Pipette 50-300 µL of the diluted HRP conjugate to each well of the ELISA at the appropriate step in the assay.
6. Store remaining conjugate solution at 2-8°C. Protect from direct exposure to light. This can be accomplished by covering the container with aluminum foil.

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.*



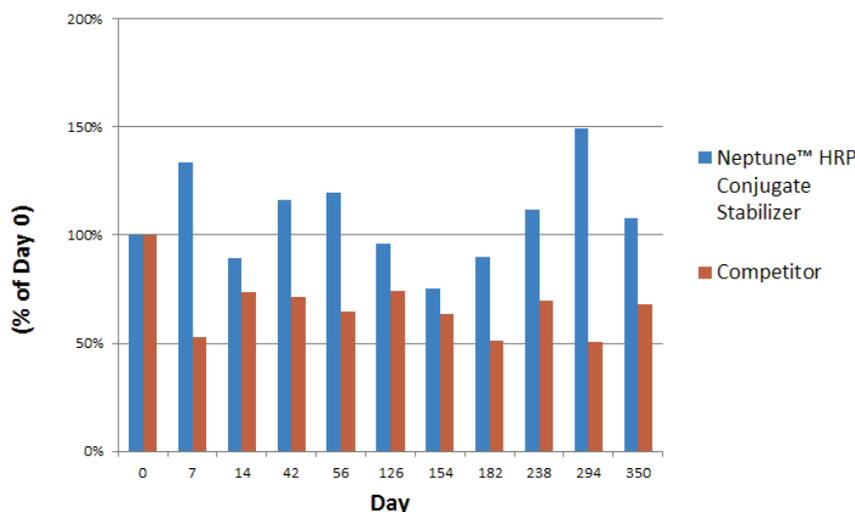
9401 James Avenue S., #155, Bloomington, MN 55431 USA  
 TOLL-FREE 1-800-829-3194 LOCAL 952-888-8788  
 FAX 952-888-8988 [WWW.IMMUNOCHEMISTRY.COM](http://WWW.IMMUNOCHEMISTRY.COM)

## Neptune™ HRP Conjugate Stabilizer, Non-Mammalian, 1X

continued from page 1...

stabilizers offers additional utility in the form of improved protein stabilization efficiency. In addition, the heterogeneous, smaller molecular weight protein matrix allows for creation of more intimate associations between stabilizer proteins and the performance-critical regions of the HRP-IgG conjugate while also serving to inhibit non-specific conjugate binding interactions with the ELISA plate surfaces. By reducing non-specific conjugate bridging events on the plate surface, undesirable background signal-related losses in sensitivity can be all but eliminated.

### SIGNAL RETENTION OF HRP-IgG CONJUGATE AT 2-8°C WITH NEPTUNE™ HRP CONJUGATE STABILIZER, NON-MAMMALIAN, 1X VS. COMPETITOR



HRP-IgG conjugate was diluted to working strength (1:60,000) in Neptune™ HRP Conjugate Stabilizer, Non-Mammalian, 1X or in a competing product and stored at 2-8°C. Performance was monitored over 12 months in an antigen-down ELISA format. Neptune™ HRP Conjugate Stabilizer, Non-Mammalian, 1X is effective at retaining both the binding and enzymatic activity of HRP conjugates. It enabled complete retention of the original signal while the signal from conjugate stored in a competing product decreased by more than 30% over the same period.

*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)