

## Signal Guard™ HRP Conjugate Stabilizer

### Ordering Information

Product Number: 11010 (50 mL)

### Storage Conditions

Store at 2 to 8 °C  
Expiration date is one year upon receipt.

### Introduction

This HRP stabilizer solution contains Tris buffered saline and proprietary reagents to stabilize and preserve Horseradish Peroxidase (HRP) conjugated primary or secondary antibodies in solution. When used with HRP conjugated primary or secondary antibodies or other proteins or compounds in solution, the HRP Conjugate Stabilizer stabilizes both the antibody and enzyme portions of the conjugate. The stabilizer prevents the loss of catalytic activity and maintains the active conformation of the antibody portion of the conjugate. With Signal Guard™ Conjugate Stabilizer, conjugates can be stored at a ready-to-use concentration without any further dilution, which avoid the risk of dilution variability for the end users. Our HRP conjugate stabilizer makes repeated assays more reproducible.

### Guidelines for Use

*Note 1: This protocol only provides a guideline, should be modified according to your specific needs.*

*Note 2: Use buffers, such as Hepes, Tris, Tricine, Mops, PIPES or MES, that contain < 10% phosphate with this HRP Conjugate Stabilizer. Buffers containing > 10% phosphate or significant amounts of borate and carbonate may cause precipitation.*

- Dilute the primary or secondary peroxidase conjugated antibody or other protein or compound with the HRP Conjugate Stabilizer to the desired 1X working concentration determined for your conjugate.
- This HRP conjugate stabilizer dilutes and stabilizes HRP conjugates in a single step.
- Do not freeze the HRP conjugates once the HRP conjugate stabilizer is added. Many HRP-conjugates prepared in this manner maintain HRP activity for up to 12 months or longer when properly stored.

### Recommended Dilution

This product is a "ready-to-use" 1X diluent for HRP conjugated primary or secondary antibody in solution. No further dilution is needed.

**Disclaimer:** *This product is for research use only and is not intended for therapeutic or diagnostic applications. Please contact our technical service representative for more information.*