

# PermaShield Clean Peroxidase Conjugate Stabilizer (RB-SPRPX)

---

## **Product Description:**

Preserves and stabilizes the activity of HRP conjugates, enzymes and other bio-molecules used in immunoassays and molecular assays. PermaShield Clean Peroxidase Conjugate Stabilizer does not contain mammalian proteins, bovine proteins, or BSA. This eliminates or minimizes background issues, unwanted protein interactions, and concerns regarding any possible transmission of bovine spongiform encephalopathy (BSE). PermaShield Clean Peroxidase Conjugate Stabilizer is not adversely affected by freezing. It provides superior HRP protection compared to competitor products. The stabilizer is supplied as a one component liquid ready to use formula.

## **Hazard Identification:**

See the material safety data sheet (MSDS).

## **Product Stability, Storage and Specifications:**

PermaShield Clean Peroxidase Conjugate Stabilizer is stable for a minimum of 2 years at room temperature. Recommended storage is at 2C to 8C. Super Shield™ can be maintained at room temperatures for extended time periods during use. Freezing will not adversely affect performance.

## **Product Use:**

PermaShield Clean Peroxidase Conjugate Stabilizer is designed as a ready-to-use buffer and stabilizer for immunoassays and molecular assays utilizing HRP labeled molecules. It can be used for high concentration stock and dilute conjugate solutions. PermaShield Clean Peroxidase Conjugate Stabilizer can be used directly as an assay buffer for incubating HRP-conjugates. It can be used for ELISA, western blots, southern/northern blots, and with lateral flow strips. PermaShield Clean Peroxidase Conjugate Stabilizer is stable and can be used at room temperature or 2C to 8C. It stabilizes dilute conjugates for months at room temperature.

These products are for research and manufacturing use only and are not intended for use in humans, therapeutic or diagnostic purposes. Sales are without any seller's warranty or representation, expressed or implied, by usage or otherwise; no claims beyond replacement of unacceptable material and no refund of purchase price shall be allowed. All claims must be made within 30 days following date of delivery.