

PermaShield Clean Universal Protein Stabilizer (RB-SPRUNV)

Product Description:

PermaShield Clean Universal Protein Stabilizer is a superior all-purpose protein, biomolecule and analyte stabilizer / diluent. It preserves and stabilizes the activity of various biomolecules used in immunoassays and molecular assays. It can be used as a stabilizer, buffer and diluent. PermaShield Clean Universal Protein Stabilizer does not contain any mammalian proteins. This eliminates or minimizes background issues, unwanted protein interactions, and concerns regarding any possible transmission of bovine spongiform encephalopathy (BSE). It is compatible and works well with RayBiotech Coat- Block reagents. PermaShield Clean stabilizer is not adversely affected by freezing. It provides superior stabilization compared to competitor products. The stabilizer is supplied as a one component liquid ready to use formula. Antimicrobial agents are included to retard bacterial growth.

Hazard Identification:

See the material safety data sheet (MSDS).

Product Stability, Storage and Specifications:

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

Product Use:

RB-SPRUNV is designed as a universal ready-to-use stabilizer and buffer for immunoassays and molecular assays. It is specially formulated to stabilize the activity of various Proteins and Biomolecules. RB-SPRUNV can be used for high concentration stock and dilute solutions. RB-SPRUNV can be used directly as an assay buffer. It can be used for ELISA, Western Blots, Southern/Northern Blots, and in lateral flow assays. It stabilizes dilute biomolecules 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.