

Secure Shield Ultra Clean Non-Milk Blocking and Diluent Solution (RB-SCRSLT)

Product Description:

Secure Shield Ultra Clean blocks and stabilizes coated ELISA plates and other solid phase substrates. RB-SCRSLT is a superior reagent that blocks, seals, and stabilizes with a consistent coat. This Coat Block was developed to substantially reduce non specific antibody binding and eliminate non-specific protein interactions. Secure Shield Ultra Clean contains no Bovine serum proteins. It provides structural support and crowding agents and minimizes background interactions and issues. RB-SCRSLT eliminates BSE regulatory concerns. RayBiotech RB-SCRSLT is compatible and works well in combination with RayBiotech Stabilizers. It is a room temperature stable reagent and stable through freeze cycles.

Hazard Identification:

See the material safety data sheet (MSDS).

Product Stability, Storage and Specifications:

Secure Shield Ultra Clean Coat Block is stable for a minimum of 2 years at room temperature. Recommended storage is at 2C to 8C. Secure Shield Ultra Clean RB-SCRSLT can be maintained at room temperatures for extended time periods during use. Freezing will not adversely affect performance.

Product Use:

RB-SCRSLT can be used for antigens or antibodies. RB-SCRSLT can be used for coating and blocking. An additional blocking step may not be required, potentially enabling blocking and stabilization in one step. Coated plates dried with RB-SCRSLT have a long term shelf life, even at room temperature, when stored in a cool and dry environment. The product is supplied as a 1X, one component liquid, ready to use formula. Antimicrobial agents are included to retard bacterial growth.

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.