

# Secure Shield Ultra Clean Non-Milk Blocking and Diluent Concentrate (RB-SCRCNC)

---

## **Product Description:**

Secure Shield Concentrate RB-SCRCNC blocks and stabilizes coated ELISA plates and other solid phase substrates. RB-SCRCNC is a superior reagent that blocks, seals, and stabilizes with a consistent coat and can be used as an excellent buffered diluent for proteins and biomolecules. This Coat Block was developed to substantially reduce non specific antibody binding and eliminate non-specific protein interactions. Secure Shield Concentrate RB-SCRCNC contains no mammalian proteins. It provides structural support and crowding agents and minimizes background interactions and issues. RB-SCRCNC eliminates BSE regulatory concerns. Data shows that Raybiotech RB-SCRCNC synergizes with RB-SPRPX PermaShield Stabilizer to further improve unwanted non specific binding. It is a room temperature stable reagent and is stable through freeze cycles.

**Hazard Identification:** See the material safety data sheet (MSDS).

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

Secure Shield Concentrate RB-SCRCNC is stable for a minimum of 2 years. Recommended storage is at 2°C to 8°C. Secure Shield Concentrate can be maintained at room temperature for extended time periods during use or when stored. Allow Secure Shield Concentrate to equilibrate to room temperature before preparing a working dilution. Freezing will not adversely affect performance.

## **Product Use:**

RB-SCRCNC can be used for antigens or antibodies. RB-SCRCNC can be used for coating, blocking and as a diluent. An additional blocking step may not be required, potentially enabling blocking and stabilization in one step. The product is supplied as a (5X-10X) concentrate. Equilibrate RB-SCRCNC to ambient temperature and dilute with RODI water to make a working solution. When diluted 5X, it serves as an excellent coat block equivalent to Raybiotech 1X Non Mammalian Protein Coat Block (RB-SCRCLN). When diluted 10X, it serves as an excellent general neutral buffered diluent. Coated plates dried with working RB-SCRCNC have a long term shelf life, even at room temperature, when stored in a cool and dry environment. Antimicrobial agents are included to retard bacterial growth.

<p>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.</p>
--