

Yttrium Silicate (YSi) SPA Reagents, Type 1 for homogeneous RIA

Product Number: RPN141 (Anti-Mouse)

Warning

For research use only.

Not recommended or intended for diagnosis of disease in humans or animals.

Do not use internally or externally in humans or animals.

Storage

Store at 2–8°C.

Expiration

Once Reconstituted, the beads are stable for up to 7 days when stored in the appropriate conditions.

Safety Warnings and Precautions

All chemicals should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.

CAUTION: For use with radioactive material.

This product is to be used with radioactive material. Please follow the manufacturer's instructions relating to the handling, use, storage, and disposal of such material.

DESCRIPTION

These generic products have been designed to enable researchers to convert existing heterogeneous assays to homogeneous scintillation proximity assay systems. A range of generic reagents, consisting

of second antibodies or protein A coupled to fluomicrospheres is provided lyophilized. Upon reconstitution, 100 µl/tube scintillation proximity assay reagent is optimal for assay performance in most assays.

REAGENT PREPARATION

Reconstitute the bottle containing scintillation proximity assay reagent with 50 ml assay buffer. Gently shake the bottle for 5 minutes to dissolve the buffer salts.

Note: The scintillation proximity assay fluomicrospheres are insoluble. The bottle contents should be magnetically stirred to ensure a homogeneous suspension when pipetting into assay tubes/wells.

Storage: After reconstitution the scintillation proximity assay reagent will be stable for up to 7 days stored at 2-8°C. For extended storage, store at -15°C to -30°C.

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

RPN141-R Rev01

The logo for Revvity, featuring the word "revvity" in a lowercase, sans-serif font.