

Phosphodiesterase SPA Beads

Product Number: RPNQ0150 (500mg containing 18mM zinc sulfate)

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 4°C to 8°C. Do not freeze.

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.

Quality Control

Each batch of PDE SPA beads is tested for functional performance using the a [3H] SPA enzyme assay.

Critical Parameters

The following points are critical

- Reconstitute to 20 mg/ml in water. Using a different volume will alter the concentration of the bead/zinc sulphate suspension and affect assay performance.
- Ensure that the beads are uniformly suspended before use by inverting the vial several times and/or placing on a vortex mixer for a few seconds.
- The beads will settle completely after standing for 15 minutes. During dispensing steps, it is recommended that the beads are stirred continuously using a magnetic stirrer to ensure an even suspension.

Counting

1. SPA PDE beads in general will yield approximately 65-70% of the cpm expected from conventional liquid scintillation counting.
2. Scintillant should not be added to the assay.
3. If assay tubes are used and the scintillation counter has removable racks, tubes can be loaded directly into the racks.
Alternatively, assay tubes can be placed into standard liquid scintillation vials and loaded into the counter.
4. For counters fitted with spectrum analysis packages the suitable window opening should be determined and the machine set accordingly. Windows for other counters should be set wide open.
5. When researchers are using highly colored samples, color quench correction may be necessary.

Please note: The SPA counts obtained will depend on the type of counter used and the absolute efficiency of the instrument.

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.

RPNQ0150-R Rev01

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a modern, clean appearance.