

Wheatgerm Agglutinin PEI Type A PS SPA Beads

Product Number: RPNQ0286 (50mg)

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

Wheatgerm agglutinin SPA imaging beads are supplied as 50 mg lyophilized solid containing 10% sucrose by weight. This material should be stored, protected from light, 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.

Quality Control

Each batch of wheatgerm agglutinin (WGA) - PEI Type A or B SPA Imaging beads is tested for its relative binding capacity of [³H] N,N',N''-triacetylchitotriose and [³H methyl,1,2] thymidine triphosphate.

BEAD RECONSTITUTION

Before use the beads should be reconstituted in a buffer appropriate for the particular assay to be performed. Standard buffers such as Tris, PBS and HEPES have all been routinely used in receptor SPA Imaging applications. The WGA-PEI Type A or B SPA Imaging beads are stable in a range of pH and solvents when used at typical assay concentrations i.e. <10% (v/v). However, it remains the responsibility of the user to evaluate the effect of the buffer composition and pH on the assay. The WGA-PEI Type A or B SPA Imaging beads should be mixed to ensure a

homogeneous suspension whilst pipetting. This may be done by either recapping and shaking the bottle at regular intervals or by gently stirring with a magnetic stirrer.

Reconstituted beads can usually be stored at 2-8°C for up to seven days. DO NOT FREEZE.

PLEASE NOTE: The WGA-PEI Type A or B SPA Imaging beads have been freeze-dried from a 1% sucrose solution. Antimicrobial agents are not included in this reagent. The user should therefore be aware that microbial contamination might occur when the reconstituted beads are stored for prolonged periods. If antimicrobial agents (e.g. Sodium Azide) are added on storage, then it remains the responsibility of the user to evaluate the effects of the added agent on the assay.

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.

RPNQ0286-R Rev01

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