

Polystyrene His Tag Imaging Beads

Product Number: RPNQ0266 (500mg)

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

PS his tag imaging beads are supplied by as an aqueous suspension at a concentration of 100 mg/ml. 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 polystyrene (PS) his tag imaging beads* is tested for its relative binding capacity of the peptide [3H]Tyr His His His-His-His-His-Ala.

BEAD RECONSTITUTION

Before use, the PS his tag imaging beads should be diluted in a buffer appropriate for the particular assay to be performed. The beads should be mixed to ensure a homogeneous suspension while pipetting. This may be done by continuous agitation with a magnetic stirrer.

Note: Magnetic stirrer bars should be coated with a chemically inert material and be free from any surface bound metals or metal salts.

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

PLEASE NOTE: Anti-microbial agents are not included in this reagent.

The user should therefore be aware that microbial contamination may occur when the diluted beads are stored for prolonged periods. If anti-microbial agents (eg 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.

ASSAY CONDITIONS

PS his tag imaging beads may be used in a direct assay format where bead is used to quantify the binding of a directly radiolabeled histidine tagged fusion protein, peptide or oligopeptide. Alternatively, PS his tag imaging beads may be used in an indirect assay format where bead is used to quantify the association of a radiolabeled binding partner to a histidine tagged fusion protein, peptide or oligopeptide. The binding of radiolabeled ligands brings the isotope into close proximity with the scintillant which is incorporated within the bead. This allows the emitted radiation (beta-particles for [3H] or Auger electrons for [125]) to stimulate the scintillant to emit light. Any unbound radiolabeled ligand is not in close enough proximity to the scintillant to allow such energy transfer and hence no signal is generated. Light emitted by stimulated LEADseeker imaging beads is detected using a Viewlux imager. Other isotopes such as [33P] and [35S] can also be used in LEADseeker imaging format. It remains the responsibility of the user to optimize the amount of reagent bead required and the incubation time required for each assay. To achieve optimal counts, excess bead should be present in order to capture all of the receptor present in the assay tube. The amount of receptor preparation together with the radiolabeled ligand being used needs to be optimized for each 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.

RPNQ0266-R Rev01

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