

GTP Agarose high (>6μmol/ml)

Applicable to: 505-0001 2ml – GTP Agarose high (>6μmol/ml)
502-0002 5ml – GTP Agarose high (>6μmol/ml)

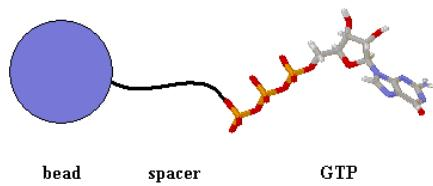
Release 1. © EXPEDEON. 28/02/2018

INTRODUCTION

Affinity resins have been widely used for the purification of enzymes and other proteins that bind nucleotides and related molecules.

GTP Agarose comprises GTP attached to agarose beads. A long hydrophilic spacer (14-atom) is used to minimise unwanted hydrophobic interactions and to facilitate unhindered interactions with biomolecules. The GTP is coupled via its -phosphate group which means that the resin is resistant to phosphatases found in many crude tissue extracts.

Figure 1. Schematic diagram of GTP Agarose



STORAGE AND SHIPPING

The resin is supplied as 50% (v/v) slurry in 10mM Tris/300mM NaCl/1mM EDTA, pH 8.0. The product is shipped at ambient temperature but should be stored at 4°C upon arrival.

MATERIALS REQUIRED (BUT NOT SUPPLIED)

For small sample volumes you may need only a microfuge and 1.5ml tubes. For larger volumes (up to 20ml) purification of binding proteins is conveniently carried out using disposable polypropylene columns. A simple mixing device (e.g. rotary shaker or end-over-end mixer) may also be useful.

OVERVIEW OF PROCEDURE

GTP Agarose is added to a crude protein extract and the suspension is gently mixed. After a period of incubation the resin is transferred to a disposable column and washed to remove non-bound or loosely adsorbed material. Finally, the column is eluted with buffer containing a competing ligand.

Since GTP Agarose may capture more than one type of GTP-binding protein the instructions below provide only general guidance on the use of resin. You may need to modify the conditions to facilitate the binding of your particular biomolecule of interest.

BUFFERS

For simplicity we would recommend that you start with the same buffer for the equilibration, binding and wash steps. The elution buffer is prepared by adding a competing ligand.

Types of buffers

The buffer and pH must be compatible with the biomolecule of interest. Tris (pH 7.5-8.5) and Hepes (pH 7.0-8.0) are commonly used but other buffers may also be suitable.

Metal ions

Binding proteins often recognise metal ion-nucleotide complexes rather than free nucleotide. It may be necessary therefore to include MgCl₂ or other suitable metal salt (usually at least 10mM) in column buffers to facilitate metal-dependent interactions with the resin.

Salts

To prevent non-specific electrostatic interactions with the matrix it is usual to include 100mM–500mM NaCl, KCl or other salt in the buffer.

Thiols

Thiols are often included in buffers to prevent oxidation of cysteine residues. A final concentration of 1mM DTT is commonly used. DTT is not stable and should be added to the buffer immediately before use.

Protease inhibitors

Protease inhibitors (e.g. PMSF, benzamidine) may or may not be required, depending on the sensitivity of the protein of interest to proteolysis. It is also advisable to carry out the binding and wash steps in a cold room or fridge using ice-cold buffers.

Detergents

Detergents (e.g. Triton X-100) are sometimes used to prevent non-specific hydrophobic interactions. Since the resin and spacer are hydrophilic a detergent may not be necessary. However, if a detergent is required try relatively low concentrations (0.02-0.1%) in the first instance.

CHROMATOGRAPHY STEPS

Make sure the resin has been fully equilibrated with the column equilibration buffer before commencing the purification procedure. Dialyse or desalt the sample into the same buffer before application to the resin.

Binding step

If you do not have access to an automated chromatography system, a batch-binding method may be used. Protein samples with volumes of 0.5-1.0ml should be incubated in 1.5ml tubes with 50-100μl of GTP Agarose. For larger sample volumes the incubation should be carried out in 10ml, 30ml or 50ml tubes (or in a capped disposable column with an integral upper reservoir). Allow at least 1 hour at 4°C for binding to take place, and agitate the sample at regular intervals to prevent settling of the resin.

If you have a pump system the recommended flow rate in the first instance is 0.1-0.25 ml/min for columns that are 1-5ml in size, though you may wish to explore higher flow rates especially if the volume of material to be processed is large.

Wash step

If incubations have been carried out in small tubes, the resin should be subjected to five or more cycles of washing and centrifugation (e.g. in a microfuge for 3-4 seconds) using ice-cold buffers. On a larger scale it is easier to transfer the suspension to a disposable polypropylene column and to allow the non-bound material to drip

through under the force of gravity. Add the wash buffer carefully down the inner surface of the column and try not to disturb the resin otherwise the wash buffer will mix with the non-bound material, leading to less efficient washing of the resin. It is important to remove all of the non-bound material prior to elution. The absence of protein in the washes is easily verified with a dye-based protein detection reagent (e.g. Bradford reagent) or with a UV monitor.

Elution step

It is important to appreciate in affinity chromatography that the eluting ligand (competing ligand or 'displacer') does not usually drive the bound protein from the resin; rather, it associates with proteins that dissociate from the resin and prevents their rebinding. The concentration of the displacer has to be sufficiently high to compete with any unoccupied ligand sites on the resin and sufficient time has to be allowed for dissociation to take place. If GTP is employed as the competing ligand a concentration in the range 5-10mM is a useful starting point.

For experiments carried out in 1.5ml tubes, the elution buffer (0.25-1.0ml) is added to the resin. After >30 minutes the resin is centrifuged and the supernatant fraction is carefully removed. If a drip column format is used, the displacer is allowed to pass into the column bed and the flow is then halted (e.g. by capping the column outlet). After a period of equilibration (15-30 minutes) the dissociated proteins are flushed out by application of more elution buffer. This step can be repeated until protein is absent from the eluted fractions. If a pump is available the column can be eluted using continuous flow at a rate of 0.05-0.1ml/min, but it may be necessary to reduce the flow rate (or switch off the pump for a period of time) to ensure that the protein elutes in a relatively small volume.

COLUMN REGENERATION

After each run, wash the column with a neutral buffer containing 1M NaCl and then re-equilibrate with 10mM Tris/30 mM NaCl/1mM EDTA pH 8.0. Do not wash the column with strong acid or base. For long-term storage add a preservative (e.g. 0.1% sodium azide).

RELATED PRODUCTS

Description	Prod. Code
Control Resin	520-0002
	520-0005

This is agarose with no GTP bound. It should be loaded, washed and eluted in exactly the same way as the GTP Agarose.

Limited use license

Expedeon's products are offered for research purposes alone, and are not intended for human, therapeutic or diagnostic use. The purchase of this product conveys to the buyer (whether the buyer is a not-for-profit, academic or for-profit entity) the non-transferable right to use the amount of product purchased and the components of the product for in-house research. The buyer shall not sell or otherwise transfer this product, its components, or materials prepared therefrom to any third party. The buyer shall not use this product or its components for commercial purposes. For the avoidance of doubt, 'commercial purposes' means any activity by a party for consideration and includes, without limitation, use of the product or its components (i) in the manufacturing of other materials (ii) to provide a service, information or data, (iii) for therapeutic, diagnostic or prophylactic purposes, or (iv) for repackaging/resale, whether or not such product or its components are resold for use in research. The use of this product by the buyer constitutes agreement with the terms of this limited use label.

Many of Expedeon's products and technologies are protected by patents or are the subject of patent filings. For information on purchasing a license for commercial use please contact us at +44 (0) 1223 873364