

Version 1a Last updated 19 August 2020

ab270533

GTP-Agarose

A product of Expedeon, an
Abcam company

Applicable to Expedeon product codes 505-0002.

View ab270533

GTP-Agarose datasheet:

www.abcam.com/ab270533

(use www.abcam.cn/ab270533 for China, or www.abcam.co.jp/ab270533 for Japan)

This product is for research use only and is not intended for
diagnostic use.

Table of Contents

1. Overview	2
2. Materials Supplied and Storage	3
3. Materials Required, Not Supplied	3
4. Technical Considerations	4
5. Chromatography Steps	6
6. Frequently Asked Questions	8

1. Overview

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

GTP-Agarose (ab270533) comprises GTP attached to agarose beads. A long hydrophilic spacer (14-atom) is used to minimize 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.

2. Materials Supplied and Storage

Store resin at +4°C. **Do not freeze. Freezing the suspension will damage the agarose beads.** The resin is supplied as 50% (v/v) slurry in 10mM Tris/300mM NaCl/1mM EDTA, pH 8.0.

Item	Quantity	Storage temperature
GTP-Agarose	2 mL	+4°C

3. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay. For small sample volumes you may need only a microfuge and 1.5 mL tubes. For larger volumes (up to 20 mL) 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.

4. Technical Considerations

4.1 Overview of the 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.

4.2 Buffer recommendations:

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 recognize metal ion-nucleotide complexes rather than free nucleotide. It may be necessary therefore to include MgCl_2 or other suitable metal salt (usually at least 10 mM) 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 100 mM–500 mM 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 1 mM 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.

5. 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.

5.1 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 0.5-1.0 mL should be incubated in 1.5 mL tubes with 50-100 μ L of GTP Agarose. For larger sample volumes the incubation should be carried out in 10 mL, 30 mL or 50 mL 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 1-5 mL in size, though you may wish to explore higher flow rates especially if the volume of material to be processed is large.

5.2 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.

5.3 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 5-10 mM is a useful starting point.
- For experiments carried out in 1.5 mL tubes, the elution buffer (0.25-1.0 mL) 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 0.05-0.1 mL/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.

5.4 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).

6. Frequently Asked Questions

6.1 What is the best buffer to use?

There is no “right” answer here. In the absence of any information on the binding requirements of the protein(s) of interest a good starting point is a buffer containing 20 mM Hepes, 100-500 mM NaCl (or KCl), 20 mM MgCl_2 and 1 mM DTT, pH 7.5. Alternatively, try experimenting with several buffer conditions using small amounts of resin in 1.5 mL tubes. After the wash step, elute with GTP or other competing ligand and analyse the eluted proteins by SDS-PAGE. The purification method can then be scaled up using the preferred buffer conditions.

6.2 Are there any other ways of eluting binding proteins from the GTP Agarose?

Yes. While GTP is the obvious choice, ligands that are structurally related to GTP may be used (e.g. GMP and guanosine) to elute a specific subset of the GTP-binding proteins. Drugs that are known to bind to GTP-binding proteins might also be used. If preservation of biological activity is not required, aliquots of resin may be boiled with SDS sample buffer prior to gel electrophoresis.

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)