

# Dynabeads™ His-Tag Isolation & Pulldown

Catalog Nos. 10103D, 10104D

Store at 2°C to 8°C

Publication No. MAN0017121

Rev. A.0

## Kit contents

Cat. No.	Volume	Capacity
10103D	2 mL	40 tests
10104D	10 mL	200 tests

Dynabeads™ His-Tag Isolation & Pulldown contains 40 mg beads/mL in 20% ethanol and has a capacity of isolating 40 µg of a 28 kDa histidine-tagged protein/ mg (25 µL) beads.

## Product description

Dynabeads™ His-tag Isolation & Pulldown is used for the isolation of histidine-tagged proteins.

The optimized cobalt-based immobilized metal affinity chromatography (IMAC) chemistry on the Dynabeads™ magnetic beads bind histidine-tagged proteins with higher selectivity than agarose- and sepharose-based bead systems.

Dynabeads™ magnetic bead-based technology makes the purification quick and easy:

Add Dynabeads™ magnetic beads to a sample containing histidine-tagged proteins and allow the proteins to bind. Isolated proteins can be left on the Dynabeads™ magnetic beads and used directly in downstream applications. Alternatively, the isolated histidine-tagged proteins can be eluted from the beads.

Elution conditions are less stringent than other technologies, yielding more functional isolated proteins. These characteristics make Dynabeads™ magnetic beads the ideal product for purifying histidine-tagged proteins expressed in *E. coli*.

## Required materials

- DynaMag™ Magnet (See [thermofisher.com/magnets](http://thermofisher.com/magnets) for recommendations on magnets appropriate for manual or automated protocols)
- Sample mixer allowing tilting and rotation of tubes (e.g. HulaMixer™ Sample Mixer)
- Test tubes and pipettes
- Buffers (see Table 1)

## Sample guidelines

- Ensure that the lysate does not contain:
  - EDTA (or other chelators)
  - Ionic detergents
  - DTT or DTE
- The lysate should have a pH between 7 and 8
- Methods for preparing cell lysates include use of:
  - 1X Binding/Wash Buffer (see “Protocol”) with 1% Triton™ X-100 (for mammalian and insect cells only)
  - French press
  - Sonication
 Efficiency of lysis can be increased by the addition of lysozyme
  - Alternative lysis strategies for *E. coli* include use of commercially available ready-made lysis buffers.

Table 1: Required buffers

2X Binding/Wash Buffer*	His Elution Buffer	2X Pull-down Buffer*	Buffer modifiers
<ul style="list-style-type: none"> <li>• 100 mM Sodium-Phosphate, pH 8.0</li> <li>• 600 mM NaCl</li> <li>• 0.02% Tween™-20</li> </ul>	<ul style="list-style-type: none"> <li>• 300 mM Imidazole</li> <li>• 50 mM Sodium-phosphate pH 8.0</li> <li>• 300 mM NaCl</li> <li>• 0.01% Tween™-20</li> </ul>	<ul style="list-style-type: none"> <li>• 6.5 mM Sodium-phosphate, pH 7.4</li> <li>• 140 mM NaCl</li> <li>• 0.02% Tween™-20</li> </ul>	<ul style="list-style-type: none"> <li>• 1 M NaCl</li> <li>• 0.1 M Imidazole</li> </ul>

\* Note that the 2X Binding/Wash Buffer and the 2X Pull-down Buffer need to be diluted to 1X concentration prior to use.

Alternative binding and/or washing buffers may also be used for isolation of your specific recombinant protein.

## General guidelines

- Add DNase I to prevent formation of a sticky pellet.
- We generally recommend applying the tube to the magnet for 2 min, but the sample can be handled when the beads are visually observed to be collected at the tube wall and the liquid is clear.
- Cell types other than *E. coli* (e.g., yeast or mammalian) can also be used for isolation of expressed histidine-tagged proteins, but optimization of the purification protocol is required.
- Protocols for the purification of histidine-tagged proteins using other metal based IMAC technologies can easily be adapted for cobalt-based IMAC, with some optimization.

## Protocol

We recommend preparing your sample containing the histidine-tagged protein in a total volume of 700 µL 1X Binding/Wash Buffer.

1. Thoroughly resuspend the Dynabeads™ magnetic beads in the vial (vortex >30 sec or tilt and rotate 5 min).
2. Transfer 50 µL (2 mg) Dynabeads™ magnetic beads to a microcentrifuge tube. Place the tube on a magnet for 2 min. Aspirate and discard the supernatant. Add your sample (prepared in 1X Binding/Wash Buffer) to beads. Mix well.
3. Incubate on a roller for 5 min at room temperature (or colder if the protein is unstable at room temperature). The incubation time may be increased up to 10 min.
4. Place the tube on the magnet for 2 min, then discard the supernatant.
5. Wash the beads 4 times with 300 µL 1X Binding/Wash Buffer by placing the tube on a magnet for 2 min and discard the supernatant. Resuspend the beads thoroughly between each washing step.
6. If the protein is to be eluted, proceed to step 7.
  - To use bead/protein complexes in other applications, resuspend the bead/protein complex in a suitable volume of 1X Pull-down Buffer (or other buffer compatible with your downstream application).
  - If you wish to continue with Pull-down, continue to step 1 in “Protein pull-down” (see page 2).
7. Add 100 µL His-Elution Buffer. Incubate the suspension on a roller for 5 min at room temperature (or colder if the protein is unstable at room temperature).
8. Apply on the magnet for 2 min and transfer the supernatant containing the eluted histidine-tagged protein to a clean tube.

## Protein pull-down

1. Prepare your sample in 1X Pull-down Buffer in a total volume of up to 700  $\mu$ L.
2. Add your sample (prepared in 1X Pull-down Buffer) to the bead/protein complex from step 5 in "Protocol" (see page 1).
3. Incubate on a roller for 10 min at room temperature (or cold if the protein is unstable at room temperature). The incubation time may be increased up to 30 min.
4. Place the tube on a magnet for 2 min, then discard the supernatant.
5. Wash the beads 4 times with 300  $\mu$ L 1X Binding/Wash Buffer by placing the tube on a magnet for 2 min and discard the supernatant. Resuspend the beads thoroughly between each washing step.
6. Add 100  $\mu$ L His-Elution Buffer. Incubate the suspension on a roller for 5 min at room temperature (or cold if protein is unstable at room temperature). Collect the beads at the tube wall using a magnet and transfer the supernatant containing the eluted histidine-tagged protein and its interacting protein to a clean tube. The elution volume may be decreased to 50  $\mu$ L.

## Automated Purification Protocols

Protein purification using Dynabeads™ His-tag Isolation and Pulldown can easily be automated on a wide variety of platforms. Automation protocols are available at: [thermofisher.com](http://thermofisher.com)

## Description of Materials

Dynabeads™ His-tag Isolation and Pulldown are uniform, superparamagnetic beads, 1  $\mu$ m in diameter, coupled with highly specific IMAC chemistry. The technology is comprised of a tetradentate metal chelator in which four of cobalt's six coordination sites are occupied. The imidazole rings of histidine residues present in a polyhistidine peptide chain are able to occupy the two remaining coordination sites, resulting in protein binding.

## Related Products

Product	Cat. No.
DynaMag™-2 Magnet	12321D
HulaMixer™ Sample Mixer	15920D

**REF** on labels is the symbol for catalog number.

## Important licensing information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Manufactured by Thermo Fisher Scientific Baltics UAB, V.A. Graiciuno 8, LT-02241 Vilnius, Lithuania. Thermo Fisher Scientific Baltics UAB complies with Quality System Standards ISO 9001 and ISO 13485.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Triton is a trademark of The Dow Chemical Company. Tween is a registered trademark of Croda Americas, PLC.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

For support visit [thermofisher.com/support](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[thermofisher.com](http://thermofisher.com)

19 May 2017

**ThermoFisher**  
SCIENTIFIC