

For life science research only.
Not for use in diagnostic procedures.



cOmplete His-Tag Purification Resin

 **Version: 04**

Content Version: November 2020

Pre-charged, ready-to-use resin for small or large-scale purification of His-tagged proteins.

Cat. No. 05 893 682 001	25 ml settled resin volume
Cat. No. 05 893 801 001	200 ml settled resin volume

Store the product at +2 to +8°C.

1.	General Information	3
1.1.	Contents	3
1.2.	Storage and Stability	3
	Storage Conditions (Product)	3
1.3.	Additional Equipment and Reagent required	3
1.4.	Application	3
2.	How to Use this Product	4
2.1.	Before you Begin	4
	Sample Materials	4
	Lysate preparation	4
	General Considerations	4
	Technical specifications	6
	Volumetric and linear flow rates	6
	Safety Information	6
	Handling requirements	6
	Laboratory procedures	6
	Waste handling	6
	Working Solution	7
2.2.	Protocols	7
	Chromatography purification	7
	Column packing for FPLC applications	7
	Purification under native conditions	8
	Purification under denaturing conditions	9
	Batch purification	10
	Cleaning	10
	Purification process optimization	11
2.3.	Parameters	12
	Affinity/Binding Capacity	12
3.	Troubleshooting	13
4.	Additional Information on this Product	14
4.1.	Test Principle	14
4.2.	Quality Control	15
5.	Supplementary Information	16
5.1.	Conventions	16
5.2.	Changes to previous version	16
5.3.	Ordering Information	16
5.4.	Trademarks	17
5.5.	License Disclaimer	17
5.6.	Regulatory Disclaimer	17
5.7.	Safety Data Sheet	17
5.8.	Contact and Support	17

1. General Information

1.1. Contents

Bottle	Label	Function / Description	Catalog Number	Content
1	cOplete His-Tag Purification Resin	<ul style="list-style-type: none"> Ready-to-use resin for the purification of His-tagged proteins. Pre-charged with Ni²⁺. Supplied as a 50% suspension in a 20% ethanol storage buffer. 	05 893 682 001	1 bottle, 25 ml
			05 893 801 001	1 bottle, 200 ml

1.2. Storage and Stability

Storage Conditions (Product)

When stored at +2 to +8°C, the product is stable through the expiry date printed on the label.

Vial / bottle	Label	Storage
1	cOplete His-Tag Purification Resin	Store at +2 to +8°C.

1.3. Additional Equipment and Reagent required

For purification under native conditions

- NaH₂PO₄
- NaCl
- Imidazole
- Ethanol

For purification under denaturing conditions

- NaH₂PO₄
- Tris-HCl*
- Urea

For cleaning

- Imidazole
- HCl
- Ethanol
- SDS*
- DTT*
- Guanidinium-HCl

1.4. Application

cOplete His-Tag Purification Resin can be used for:

- Small or large-scale purification of His-tagged proteins, yielding highly purified proteins from crude lysates.
- Batch or column chromatography procedures and allows for target protein purification using both native and denaturing conditions.

i *It is compatible with common reducing agents and chelators while preventing contamination of the protein preparation with heavy metals.*

2. How to Use this Product

2.1. Before you Begin

Sample Materials

Lysate preparation

Prior to the lysate preparation, express the target protein recombinantly in a host organism, preferably in *E. coli*. The choice of an *E. coli* strain, expression conditions, such as temperature, media composition, induction strength, duration of induction, and the choice of lysis buffer can have a significant impact on the target protein yield and purity. These parameters must be optimized on a case-to-case basis.

i *Optimal methods for lysate preparation may significantly differ between different host organisms.*

⚠ *After harvesting, handle cells and lysates on ice or at +2 to +8°C.*

General Considerations

The purification of a single type of protein from a complex mixture includes a series of procedures. Consider the following variables when optimizing the protocols:

Variable	Description and Recommendation
Appropriate host organism	<p>Most proteins are produced in <i>E. coli</i> which is often the host of choice as it combines easy strain construction, rapid growth, and inexpensive culture handling with high overexpression capabilities and low background binding of host proteins.</p> <ul style="list-style-type: none"> ▪ <i>E. coli</i> features cytosolic and periplasmic expression and can be used to express recombinant proteins in both reducing and oxidizing environments. ▪ Specialized mammalian-, insect-, and yeast-based expression systems are also available, and can be used when folding problems or eukaryotic posttranslational modification problems are encountered in <i>E. coli</i>.
Native versus denaturing purifications	<p>The resin can be used to purify proteins using both native and denaturing conditions.</p> <ul style="list-style-type: none"> ▪ In most cases, a native purification is preferred as the vast majority of functional applications rely on a properly folded protein. ▪ Purification under denaturing conditions may be preferred when proteolytic degradation occurs during purification, contamination is likely from host proteins, and when low target protein solubility is present. Typical denaturing agents for such situations are urea and guanidinium-HCl. <p>i <i>The binding capacity and binding kinetics of the affinity matrix can differ in native and denaturing purification protocols.</i></p>
Appropriate length of the His-Tag	<p>The length of the His-tag fused to the target protein determines the binding specificity.</p> <ul style="list-style-type: none"> ▪ When using hexahistidine tags, histidine amino acids from the tag may not be readily distinguished from endogenous histidine amino acids. ▪ His10- to His14-tagged proteins can be more efficiently separated from histidine-rich host proteins. <p>i <i>cOmpete His-Tag Purification Resin is compatible with a polyhistidine tag length from 6 to 14 histidines.</i></p>

Appropriate buffer system	<p>The resin is compatible with a wide range of buffers. To maximize purification effectiveness, it is important to select the optimal buffer for the stability and solubility of the target protein.</p> <ul style="list-style-type: none"> Due to the high binding strength of Ni²⁺ to cOComplete His-Tag Purification Resin, the optimal buffer can be selected for protein purification without having to compromise between protein stability and resin stability. The buffer composition can also be adjusted according to the needs of the target protein. Buffers containing EDTA and DTT are compatible with cOComplete His-Tag Purification Resin. These features effectively inhibit metalloproteases and facilitate the purification of proteins prone to oxidation. To achieve optimal protein purity, the stringency during the binding and washing steps, as well as the conditions for elution, can be fine-tuned by adjusting either the concentration of imidazole or the pH value. Imidazole competes with His-tagged proteins to bind Ni²⁺ immobilized on the cOComplete His-Tag Purification Resin. Adding low concentrations of imidazole may help to revert undesired binding of host proteins to the resin. The binding of the His-tagged target protein to cOComplete His-Tag Purification Resin is also pH dependent. <p><i>i</i> The pK_a value of commonly used buffers changes with the temperature. Adjust the pH value of buffers at the same temperature as the temperature of the purification experiment.</p> <p><i>i</i> <i>E. coli</i> grown with a fermentable carbon source, such as glucose may produce organic acids that lower the pH value. Buffer the media with, for example, K_2HPO_4 and resuspend cells in an alkaline buffer with high buffering capacity. For best results, control the pH value after cell resuspension and readjust if necessary.</p>
Imidazole concentration for load/wash	<p>Nonspecific binding of proteins without a His-tag to cOComplete His-Tag Purification Resin is low. Use up to 5 mM imidazole in load and/or wash buffers.</p> <p>⚠ Do not use imidazole for new purification assays of His-tag proteins using cOComplete His-Tag Purification Resin.</p> <p>To improve the purity of the His-tag protein following this first step, use imidazole in a final concentration of up to 5 mM in a second step.</p>
Imidazole concentration for elution	<p>Up to 500 mM</p> <p><i>i</i> In contrast to other available resins, bound His-tagged protein typically elutes from the resin with a lower imidazole concentration, such as 25 to 45 mM.</p>
Compatibility for long-term storage	20% ethanol, pH 4.0 to pH 9.0
Compatibility during chromatography	The resin is compatible with 10 mM EDTA, 10 mM DTT during the purification (1 hour incubation), 6 M guanidinium-HCl, 8 M urea, pH 2.0 to 14.0.
Compatibility during cleaning	4% SDS

2. How to Use this Product

Technical specifications

Specification	Value
Matrix	Sepharose-CL 6B
Bead size	45 – 165 µM
Maximal linear flow rate	1,420 cm/hour
Recommended volumetric flow rate	The volumetric flow rate is a function of the cross section of the column. i Using the following formula, a linear flow rate can be converted to a volumetric flow (ml/min): <i>Linear flow rate (cm/hour) × column cross sectional area (cm²)/60.</i> i The column cross sectional area is defined as $\pi \times r^2$, whereas π is the constant pi, and r is the inner radius of the column.

Volumetric and linear flow rates

The following table shows the flow rates for a column with a diameter of 0.64 cm.

Volumetric flow rate [ml/min]	Linear flow rate [cm/hour]
0.9	168
1.8	336
2.4	448
3.2	598
4.0	747
4.9	915
5.8	1,084
6.7	1,252
7.6	1,420
8.5	1,589
9.4	1,697
10.0	1,805

Safety Information

Handling requirements

Always maintain cOmplete His-Tag Purification Resin in buffer; do not allow it to dry.

Laboratory procedures

- Handle all samples as if potentially infectious, using safe laboratory procedures. As the sensitivity and titer of potential pathogens in the sample material varies, the operator must optimize pathogen inactivation by the Lysis / Binding Buffer or take appropriate measures, according to local safety regulations.
- Do not eat, drink or smoke in the laboratory work area.
- Do not pipette by mouth.
- Wear protective disposable gloves, laboratory coats and eye protection, when handling samples and kit reagents.
- Wash hands thoroughly after handling samples and reagents.

Waste handling

- Discard unused reagents and waste in accordance with country, federal, state, and local regulations.
- Safety Data Sheets (SDS) are available online on dialog.roche.com, or upon request from the local Roche office.

Working Solution

Buffer	Composition	For use in...
Buffer A	<ul style="list-style-type: none"> ▪ 50 mM NaH₂PO₄, pH 8.0 ▪ 300 mM NaCl 	<ul style="list-style-type: none"> ▪ Purification under native conditions. ▪ Purification under denaturing conditions (optional). ▪ Batch purification
Buffer B	<ul style="list-style-type: none"> ▪ 50 mM NaH₂PO₄, pH 8.0 ▪ 300 mM NaCl ▪ 250 mM imidazole 	<ul style="list-style-type: none"> ▪ Purification under native conditions. ▪ Batch purification
Buffer C	<ul style="list-style-type: none"> ▪ 100 mM NaH₂PO₄ ▪ 10 mM Tris-HCl* ▪ 8 M urea ▪ pH 8.0 	<ul style="list-style-type: none"> ▪ Purification under denaturing conditions. ▪ Batch purification
Buffer D	<ul style="list-style-type: none"> ▪ 100 mM NaH₂PO₄ ▪ 10 mM Tris-HCl* ▪ 8 M urea ▪ pH 6.3 	Purification under denaturing conditions.
Buffer E	<ul style="list-style-type: none"> ▪ 100 mM NaH₂PO₄ ▪ 10 mM Tris-HCl* ▪ 8 M urea ▪ pH 5.9 	Purification under denaturing conditions.
Buffer F	<ul style="list-style-type: none"> ▪ 100 mM NaH₂PO₄ ▪ 10 mM Tris-HCl* ▪ 8 M urea ▪ pH 4.5 	Purification under denaturing conditions.

2.2. Protocols

Chromatography purification

Purifications with cOComplete His-Tag Purification Resin are compatible with various standard purification formats, including batch processing, column purification with gravity flow, and automated processing using an automated chromatography system.

Column packing for FPLC applications

- 1 Resuspend cOComplete His-Tag Purification Resin by inverting the bottle several times.
- 2 Pipette the appropriate amount of resin to a chromatography column.
⚠ **Ensure that no air bubbles get trapped in the adapter or gel bed.**
- 3 Allow for the resin to settle.
- 4 Drain the excess buffer through the column by gravity flow.
- 5 Insert the top adapter and adjust it to the top of the bed.
- 6 Connect the column to a chromatography system/buffer reservoir.
- 7 Use a flow rate as described in section, **General Considerations, Technical Specifications.**

2. How to Use this Product

Purification under native conditions

The purification of native proteins requires optimal buffer conditions for the target protein. Well established buffers which can be adapted to achieve optimal conditions for a specific target protein are described. cOmplete His-Tag Purification Resin offers full flexibility in selecting the optimal buffer conditions without compromises.

⚠ cOmplete His-Tag Purification Resin has been optimized using Buffer A and Buffer B. Other buffers may function as well, but need to be tested prior to use.

i See section, **Working Solution** for additional information on preparing solutions.

1 Equilibrate the column with 10 column volumes of Buffer A.

2 Load the cleared sample containing the His-tagged protein, for example, after an ultracentrifugation or filtration step, onto the column with a volumetric flow rate of 2.5 ml/min for 5 ml bed volume of resin or 0.5 to 1 ml/min for 1 ml bed volume.

⚠ To prevent blockage of the resin, remove insoluble material prior to loading the column.

⚠ Since the binding specificity of the resin is high, the kinetics of adhesion of the protein to the resin is slower than other available resins. If using high volumetric flow rates for loading, protein yield can decrease.

3 Wash the column with Buffer A until the absorption (A_{280}) reaches the baseline level, approximately 10 column volumes.

4 Elute the His-tagged protein with a gradient of Buffer A (without imidazole) and Buffer B (250 mM imidazole).

⚠ Protein peaks can be expected between 25 to 45 mM imidazole. Due to the specific characteristics of cOmplete His-Tag Purification Resin, a protein can already be eluted with approximately 25 mM imidazole.

⚠ The amount of imidazole required for efficient release of the target protein depends on the length and accessibility of the His-tag.

5 Wash and equilibrate for the next run, see section, **Cleaning Protocols**.

⚠ If the purification column is not immediately reused, clean the column with 2 column volumes of 2 M imidazole to remove nonspecific binding of proteins. Equilibrate the column in a 20% ethanol solution, and store at +2 to +8°C to prevent cell growth.

i See sections, **Purification process optimization** and **Troubleshooting** for additional information on optimizing the purification results.

Purification under denaturing conditions

Denature the protein or dissolve the inclusion bodies in a buffer containing 6 M guanidinium-HCl or 8 M urea.

⚠ *The addition of urea to buffered solutions will cause the pH to drop. It is essential to adjust the pH of the buffer with NaOH after urea addition.*

⚠ *The binding capacity may also drop significantly if the buffer composition is suboptimal.*

i *For best results, perform an overnight incubation to bind denatured target proteins more efficiently to the resin.*

⚠ *cOmplete His-Tag Purification Resin has been tested with Buffers C, D, E, and F. Other buffers may function as well, but need to be tested prior to use.*

i *See section, Working Solution for additional information on preparing solutions.*

1 Equilibrate the column with 10 column volumes of Buffer C.

2 Load the cleared sample containing the His-tagged protein, for example, after an ultracentrifugation or filtration step, onto the column with a volumetric flow rate of 0.5 to 1 ml/min for 1 ml bed volume of resin.

⚠ *To prevent blockage of the resin, remove insoluble material prior to loading the column.*

⚠ *Since the binding specificity of the resin is high, the kinetics of adhesion of the protein to the resin is slower than other available resins. If using high volumetric flow rates for loading, protein yield can decrease.*

3 Wash the column with Buffer C until the absorption (A_{280}) reaches the baseline level, approximately 10 to 20 column volumes.

4 Wash with 10 to 20 column volumes of Buffer D.

5 Wash with 10 to 20 column volumes of Buffer E.

6 Elute with 10 to 20 column volumes of Buffer F.

7 Wash and equilibrate for next run under denaturing conditions with Buffer C or wash with Buffer A to remove the denaturing agents if the column will next be used under native conditions.

⚠ *If the column is not immediately reused, clean the column with 2 column volumes of 2 M imidazole. Equilibrate the column in a 20% ethanol solution and store at +2 to +8°C to prevent cell growth.*

i *The elution can also be performed with a gradient up to 250 mM imidazole solution instead of the pH shift option.*

i *See sections, Purification process optimization and Troubleshooting for additional information on optimizing the purification results.*

2. How to Use this Product

Batch purification

The batch purification protocol can be conducted under native conditions as well as under denaturing conditions.

- 1 Transfer the appropriate amount of cOComplete His-Tag Purification Resin to a graduated container or an empty column.

- 2 Equilibrate the column with the resin with approximately 20 column volumes of Buffer A.

- 3 Add an appropriate amount of cleared lysate to the prepared resin slurry.
 - Swirl gently and incubate the mixture at +2 to +8°C for 2 to 12 hours on a shaker.

- 4 Load the resin-lysate mixture onto an appropriate sized column.

- 5 Collect the column flow through.

- 6 Wash the column with at least 5 column volumes of Buffer A.

- 7 Elute the protein with at least 5 column volumes of Buffer B.

- 8 Wash and equilibrate with Buffer A for the next run.

⚠ If the column is not immediately reused, equilibrate the column in a 20% ethanol solution and store at +2 to +8°C to prevent cell growth.

i See sections, **Purification process optimization** and **Troubleshooting** for additional information on optimizing the purification results.

Cleaning

cOComplete His-Tag Purification Resin can be used multiple times without loss of binding capacity. Over time, some protein aggregates might accumulate leading to a decrease in efficiency of the resin. This is identified by a slower flow rate or a higher back pressure.

The cleaning procedures remove aggregates for further use of the resin. Different cleaning procedures can be carried out, based on the different applications of the resin. Once the cleaning procedure is completed, transfer the resin to 20% ethanol.

Stringent native cleaning

This method is used when non-aggregating proteins have been purified and if the column is used again for purifying the same protein.

- 1 Wash with 10 column volumes of 1 M imidazole/HCl, pH 7.5.

- 2 Wash with 10 column volumes of 4 M imidazole/HCl, pH 7.5.

- 3 Equilibrate the column with binding buffer and proceed to the next round of purification or transfer the material to 20% ethanol.

Denaturing cleaning with SDS

This method is used to remove aggregated proteins and lipids.

⚠ Perform this cleaning at +15 to +25°C to achieve the optimal solubility of SDS.

⚠ Avoid using K⁺ in this buffer to prevent precipitation with SDS.

i The SDS buffer may also contain 50 mM DTT*.

- 1 Wash with 10 column volumes of 1 M imidazole/HCl, pH 7.5.

- 2 Wash twice with 10 column volumes of 1 M imidazole/HCl, pH 7.5, 20% ethanol, 2 to 4% SDS*.

- 3 Remove SDS with 3 times 10 column volumes of 20% ethanol.

Denaturing cleaning with guanidinium-HCl

This method is used to remove aggregated proteins.

i The guanidinium-HCl buffer may also contain 50 mM DTT.

- 1 Wash with 10 column volumes of 1 M imidazole/HCl, pH 7.5.

- 2 Wash 2 times with 10 column volumes of 6 M guanidinium-HCl, 1 M imidazole, pH 7.5.

- 3 Wash 2 times with 10 column volumes of 20% ethanol.

i The choice of cleaning method depends on the protein type.

i The denaturing cleaning procedure with guanidinium-HCl presents fewer constraints than the denaturing cleaning method with SDS.

Purification process optimization

The parameters allowing for the maximal protein yield and purity might vary significantly depending on the characteristics of a given target protein. For best results, optimize the key parameters in small-scale trial purifications. To optimize the protein purification procedure for highest protein purity, determine the optimal operating conditions of the resin for the specific target protein.

- Both purity and yield of a protein preparation depends on the amount of cComplete His-Tag Purification Resin used for binding. If the amount of resin is too high in relation to the amount of available target protein, the remaining binding sites on the resin may enable background binding of lysate components. If the amount of resin is too low, the binding capacity of the resin may not be sufficient to bind all target protein, resulting in a suboptimal protein yield.
- Optimal results are obtained when the capacity of the resin matches the amount of target protein.

The capacity for a given target protein depends on several factors such as target protein size, conformation, and multimerization status, length and accessibility of the His-tag, expression level, and solubility of the His-tagged protein, lysate concentration, as well as the buffer pH and composition. For best results, determine the optimal volume of resin required for the purification of a specific protein of interest. Based on the expression rate of the protein and the volume of lysate available, determine the optimal volume of resin by performing the following pretrial:

- 1 Incubate a small quantity of settled resin with varying volumes of lysates, in parallel test experiments.

- 2 Wash the resin and elute the bound proteins.

- 3 Determine the amount of target protein in the unbound fractions and in the eluate by SDS-PAGE.

2. How to Use this Product

- 4 The volume of lysate/volume of resin ratio is optimal when only a small amount of target protein remains in the lysate and the maximal amount of protein is detected in the eluate fractions.
- For example, for the purification of 20 to 40 mg T7 RNA polymerase (97 kDa), the amount of resin needed is 1 ml.

i The yield of the target protein can be optimized by allowing more time for the protein to bind to the resin. This can be performed by reducing the flow rate during the loading step of the chromatography purification. Alternatively, batch adsorption can be carried out for up to 12 hours during a batch purification procedure, without impacting the protein integrity. For example, the adsorption of the His₆-tagged T4 gene 32 protein has been measured during a batch purification procedure after different incubation times of the protein solution with cComplete His-Tag Purification Resin in a tube, on a roller, see Figure 1.

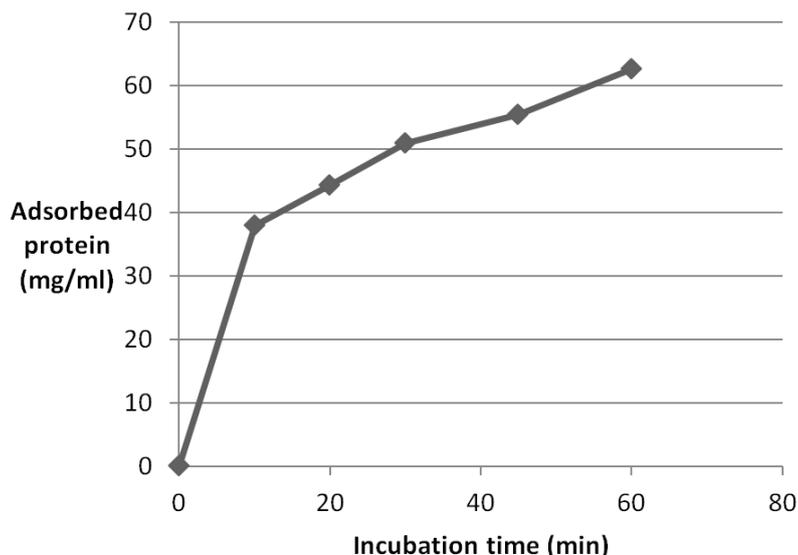


Fig. 1: Kinetics of a protein adsorption during a batch purification procedure.

- i* The optimal concentration of imidazole during binding, washing, and elution steps can be determined during pretrial experiments with a fixed aliquot of resin. This identifies the optimal amount of lysate for saturating using varying concentrations of imidazole.
- i* Optimal results can typically be achieved with buffers containing a high salt concentration (300 mM) at pH 8.0 for target proteins compatible with those conditions.

2.3. Parameters

Affinity/Binding Capacity

≥40 mg protein per ml.

- i* The binding capacity of the resin to various types of proteins may vary according to the protein characteristics, such as the size of the protein.

cComplete His-Tag Purification Resin binds with a high specificity to the polyhistidine-tagged protein.

- As a consequence, the binding kinetics may appear to be different when compared to conventional metal chelate matrices.
- Full capacity of the resin can be achieved by allowing more time for the protein to bind to the resin by lowering the flow rate during the chromatography purification procedure or by increasing the incubation time during the batch purification procedure.

3. Troubleshooting

Observation	Possible cause	Recommendation
Bubbles form in the bed resin.	Mixing of the storage buffer (20% ethanol) with aqueous buffer.	After storage at +2 to +8°C, equilibrate the resin to +15 to +25°C prior to packing the column. Degas the buffer prior to equilibration of the column.
The sample does not flow easily through the resin (low flow rate or high back pressure).	Particulates from the lysates may have clogged the resin.	Centrifuge (ultracentrifuge) the sample prior to loading on the column. Reduce the flow rate. Clean the resin using a denaturing cleaning procedure.
Inefficient binding of the target protein to the resin.	Suboptimal buffer conditions during the binding step.	Lower the imidazole concentration and/or increase the pH during the binding step.
	Incubation time is too short.	Extend the incubation time. Lower the flow rate during binding.
	The His-tag is not accessible.	Change the position of the His-tag. Use a longer His-tag.
Inefficient or no elution of the target protein.	The target protein multimerizes and binds more avidly to the resin.	Increase the imidazole concentration during elution.
	The protein precipitates on the resin prior to elution.	Increase ionic strength to minimize isoelectric precipitations. Elute under denaturing conditions.
	The target protein precipitates during a pH shift elution.	Elute with imidazole instead.
Recovery of the target protein is too low.	The target protein may be degraded.	Add protease inhibitors to the sample if degradation occurs during cell lysis. Work at +2 to +8°C.
	The His-tag might not be accessible.	Use a longer His-tag. Check if the target protein contains the His-tag. Optimize expression conditions and buffers. Change the localization of the His-tag.
	The His-tag might have been digested by proteases.	Change to another expression host. Use protease inhibitors.
	The target protein might not be soluble.	Lower the expression temperature, strength, and duration of induction. Purify under denaturing conditions. Include solubility-enhancing fusion partners.
	The resin is limiting.	Verify that the resin bed volume is proportionate to the level of expressed His-tagged protein.
	Target protein elutes with contaminants.	The host proteins interact with the resin.
DNA and/or RNA contaminants.		Wash the column with a stringent buffer. Purify under denaturing conditions. Include a DNase I digestion step and/or a Polymin P-mediated precipitation step prior to adding the lysate to the resin.

4. Additional Information on this Product

Target protein is degraded during or following the cell lysis.	Insufficient protection from proteases.	Add protease inhibitors to the buffers and/or culture.
		Optimize the experimental workflow.
		Strictly work on ice.
Target protein is degraded in the host cell.	Wrong host strain.	Use a protease-deficient host strain.
	Induction time too long.	Reduce the induction time.

4. Additional Information on this Product

4.1. Test Principle

Recombinant protein expression

Purifying a protein of interest is often essential for determining its function, structure, or interactions, for raising specific antibodies, or preparing enzymes for practical applications. Isolation of naturally expressed proteins from their original source can be a complex process involving numerous chromatographic steps. Recombinant protein expression in dedicated host organisms can greatly simplify this task. Such expression systems generally ensure higher expression levels. Fusing the target protein to a tag also confers advantageous binding ability to an affinity matrix.

Protein purification using immobilized Ni²⁺

The most common technique for efficiently obtaining large yields of highly purified protein in a short timeframe involves engineering a polyhistidine tag into the protein of interest, followed by purification using Immobilized Metal Ion Affinity Chromatography (IMAC).

- The most commonly used tag for large amounts of highly purified protein is a poly-histidine tag (His-tag). This tag has 6 to 14 histidines, typically fused to the N- or C-terminal end of a target protein.
- In some cases, the tag is also inserted into an exposed loop of the target protein.
- The imidazole side chains of a His-tag can form reversible coordinative bonds to divalent metal ions, such as Ni²⁺, Co²⁺, or Zn²⁺. This property can be used to separate polyhistidine-tagged target proteins from other proteins. Ni²⁺ show the highest affinity and selectivity for His-tags, and are therefore the preferred ions. Using a specific chelator covalently linked to a matrix, Ni²⁺ are immobilized to still permit interactions with histidine side chains. When His-tagged proteins are applied to such a Ni²⁺ resin, they specifically bind to the resin via Ni²⁺, while most untagged proteins do not. Bound proteins are released from the resin using mild conditions. Imidazole competes for coordination sites on Ni²⁺ and therefore displaces His-tagged proteins from the resin. Alternatively, lowering the pH will protonate His-tags, decreasing their affinity for the resin and hence elute the His-tagged proteins.

His-tags

Ideally, the His-tagged target protein binds much stronger to the Ni²⁺ chelate matrix than endogenous histidine-containing protein of the expression host. Relative binding strength depends on how many histidines can bind simultaneously to the matrix (avidity effect). Longer His-tags confer stronger binding and better separation of the target from potentially contaminating host proteins. The classic His-tag has six consecutive histidines. Tags with 10 to 14 histidines may produce a better purification. Most importantly, His-tagged proteins can be purified using Ni²⁺ chelate matrices under both native and denaturing conditions. Due to their hydrophilic and flexible nature, these matrices increase the solubility of the target proteins and only rarely interfere with protein function. This unique combination of features enables the His-tag to be a versatile tool for a wide range of protein purification applications.

Properties

cComplete His-Tag Purification Resin is a sepharose-based, pre-charged, ready-to-use Ni²⁺ chelate resin for small and large-scale purification of His-tagged proteins. It allows for the production of highly pure proteins from crude lysates, using a one-step purification process.

cComplete His-Tag Purification Resin is based on a chelator chemistry enabling an extremely tight binding of Ni²⁺ to the resin. In contrast to conventional nitrilotriacetic acid (NTA)-based resins and iminodiacetic acid (IDA)-based matrices, the chelator of cComplete His-Tag Purification Resin protects Ni²⁺ effectively against reduction by thiols, resulting in minimal leaching of the ions, see Figure 2. As a result, buffers and proteins processed using cComplete His-Tag Purification Resin show almost no Ni²⁺ content that would otherwise catalyze oxidative damages of proteins. As a consequence, no removal or reloading steps of Ni²⁺ are required when using cComplete His-Tag Purification Resin.

In contrast to conventional metal chelate matrices, cComplete His-Tag Purification Resin is fully compatible with solutions containing metal-chelating protease inhibitors, such as EDTA, as well as with solutions containing reducing agents such as DTT (Figure 2).

1 ml of the 3 following resins was incubated for 1 hour in a buffer containing 10 mM EDTA, 10 mM DTT, 500 mM imidazole, 300 mM NaCl, 50 mM NaH₂PO₄, pH 8.0:

- Resin 1: Commercially available Ni-NTA based resin.
- Resin 2: Commercially available Ni-Chelator based resin.
- Resin 3: cComplete His-Tag Purification Resin.

Ni²⁺ content of the buffer and of the resins was measured by ICP-MS (Inductively-Coupled-Plasma Mass-Spectrometry) to determine the amount of Ni²⁺ lost by the resins.

cComplete His-Tag Purification Resin shows <1% of Ni²⁺ content loss. It can therefore be reused and does not require to be recharged with Ni²⁺.

Resin 2 lost more than half of its Ni²⁺ content, and Resin 1 lost more than 2/3 of its Ni²⁺ content. These two Ni²⁺ based resins lost protein binding capacity and have to be regenerated but they also contaminated the protein sample with Ni²⁺.

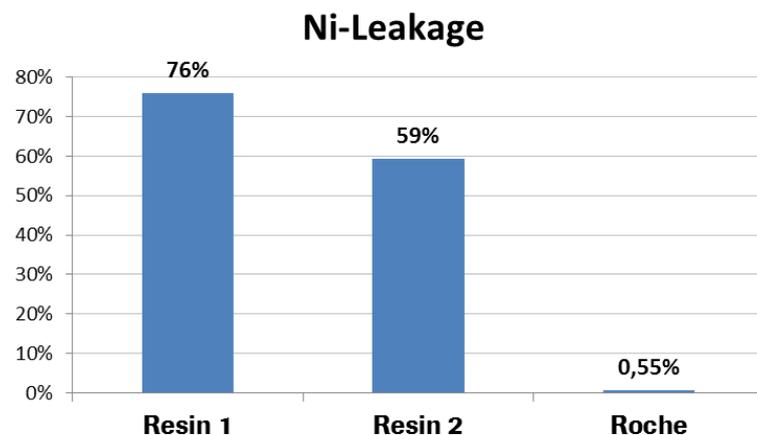


Fig. 2: Leaching of metal ions during treatment with EDTA, DTT, and imidazole.

cComplete His-Tag Purification Resin permits multiple uses (up to 5 times) without loss of binding capacity. It is compatible with solutions containing metal-chelating protease inhibitors such as EDTA, as well as with solutions containing reducing agents, such as DTT.

Together, these features enable the researchers to adapt the buffer scheme to the specific needs of the target protein over a wide range of parameters. Specifically, cComplete His-Tag Purification Resin allows for efficient protection of the target proteins from proteolytic degradation, oxidative damages, and heavy metal contamination.

The reagent compatibility also allows the resin to be used with cComplete ULTRA Tablets, Protease inhibitor cocktails containing EDTA, and with PhosSTOP Tablets for inhibiting phosphatases, as well as with common reducing agents, such as DTT.

4.2. Quality Control

For lot-specific certificates of analysis, see section **Contact and Support**.

5. Supplementary Information

5.1. Conventions

To make information consistent and easier to read, the following text conventions and symbols are used in this document to highlight important information:

Text convention and symbols

i *Information Note: Additional information about the current topic or procedure.*

⚠ Important Note: Information critical to the success of the current procedure or use of the product.

① ② ③ etc. Stages in a process that usually occur in the order listed.

① ② ③ etc. Steps in a procedure that must be performed in the order listed.

* (Asterisk) The Asterisk denotes a product available from Roche Diagnostics.

5.2. Changes to previous version

Layout changes.

Editorial changes.

Update to include new safety Information to ensure handling according controlled conditions.

5.3. Ordering Information

Product	Pack Size	Cat. No.
Reagents, kits		
Tris hydrochloride	500 g	10 812 846 001
1,4-Dithiothreitol	2 g	10 197 777 001
	10 g	10 708 984 001
	25 g	11 583 786 001
Sodium Dodecyl Sulfate (SDS)	1 kg	11 667 289 001

5.4. Trademarks

PHOSSTOP is a trademark of Roche.

All other product names and trademarks are the property of their respective owners.

5.5. License Disclaimer

For patent license limitations for individual products please refer to:

List of biochemical reagent products.

5.6. Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

5.7. Safety Data Sheet

Please follow the instructions in the Safety Data Sheet (SDS).

5.8. Contact and Support

To ask questions, solve problems, suggest enhancements or report new applications, please visit our **Online Technical Support Site.**

To call, write, fax, or email us, visit **sigma-aldrich.com**, and select your home country. Country-specific contact information will be displayed.

