

Pure Attraction

Selectivity that's simply captivating — Bio-Rad's Profinity™ IMAC Ni-charged resin provides optimal purification of recombinant His-tagged proteins.

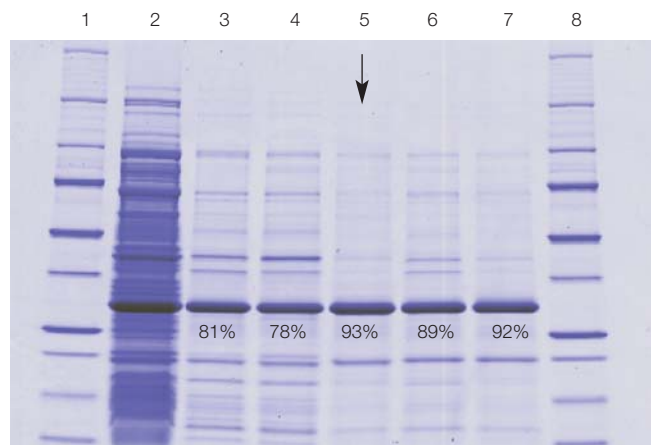
- Optimal ligand density for higher purity of target protein
- Compatibility with denaturants, detergents, and reducing agents allows excellent purification over an expanded range of conditions
- Stability in pH 1–14 for storage in a variety of solutions
- Profinity polymer bead allows purification at fast flow rates
- Easy-to-pack resin may be used in medium-pressure, gravity-flow, and spin columns
- Also available uncharged



Profinity IMAC Ni-Charged Resin

Profinity IMAC Ni-charged resin provides purification of recombinant His-tagged proteins for a wide molecular weight range. The Profinity IMAC bead is a porous 60 µm particle derivatized with iminodiacetic acid (IDA), which functions as the chelating ligand. The chemical structure of IDA, when charged with Ni²⁺ or other transition metal ions, allows highly selective binding of recombinant His-tagged proteins over naturally occurring His-containing proteins. The polymeric nature, optimized IDA ligand density, and open pore structure of the Profinity IMAC bead result in superb mechanical strength, high selectivity for target proteins, low nonspecific binding, and the ability to perform purifications at faster flow rates.

Profinity IMAC resin is stable across the full pH range (1–14) and is compatible with reagents traditionally used in the purification of His-tagged proteins. Available either charged with nickel or uncharged in small and large volumes, the resin is easy to pack in Bio-Scale™ medium-pressure, Econo-Column® gravity-flow, and Bio-Spin® columns.



Purification of a putative aminopeptidase protein using different IMAC resins. An insoluble 32 kD protein obtained from *Anabaena* sp. strain PCC 7120 (courtesy of Dr Ray Stevens, University of California, Berkeley, CA, USA) was expressed in *E. coli* and purified under denaturing conditions. *E. coli* lysate was loaded onto Micro Bio-Spin™ columns containing individual IMAC resins and purified. The binding buffer was 50 mM potassium phosphate, 300 mM NaCl, 8 M urea (pH 8.0), and the elution buffer was binding buffer plus 250 mM imidazole. To determine purity of the target protein, 3 µg of sample eluate from each column was loaded and separated on a Criterion™ gel, stained with Coomassie Blue, then quantitated using Quantity One® software. Lanes 1 and 8, 10 µl Precision Plus Protein™ standards; lane 2, 3 µl crude lysate; lane 3, Ni-charged, high-binding-capacity agarose-based resin from supplier A (IDA ligand); lane 4, uncharged agarose-based resin from supplier A (IDA ligand), charged with Ni²⁺; lane 5, Profinity IMAC Ni-charged resin; lane 6, Ni-charged agarose-based resin from supplier B (NTA ligand); lane 7, Co²⁺-charged tetradentate agarose from supplier C. The purity obtained for each resin is indicated; arrow highlights result obtained with Profinity IMAC resin.



Specifications

| | |
|-------------------------------|---|
| Functional ligand | IDA |
| Base bead | UNOsphere™ base matrix |
| Form | 50% suspension in 20% EtOH, precharged with Ni ²⁺ or uncharged |
| Particle size | 45–90 µm |
| Mean particle size | 60 µm |
| Metal ion capacity | 12–30 µmol Cu ²⁺ /ml |
| Dynamic binding capacity* | ≥15 mg/ml |
| Recommended linear flow rate | ≤600 cm/hr at 25°C |
| Maximum operating pressure | 7.5 bar (109 psi) |
| pH stability, uncharged resin | 1–14 (up to 200 hr) |
| Chemical compatibilities | See bulletin 3193 for complete list |
| Storage | 4°C to ambient temperature |
| Shelf life in 20% EtOH | >1 year at ambient temperature |
| Operational temperature | 4–40°C |
| Autoclaving conditions | 0.1 M sodium acetate at 120°C for 30 min |

* Binding capacity was determined by Q_{10%} determination under the following conditions (dynamic binding capacity will vary from protein to protein):

| | |
|----------------|--|
| Column volume | 1 ml (7 mm ID x 2.6 cm) column |
| Sample | 1.8 mg/ml pure 32 kD His-tagged protein |
| Flow rate | 1 ml/min loading 2 ml/min wash and elution |
| Loading buffer | 50 mM sodium phosphate, 300 mM NaCl, 5 mM imidazole (pH 8.0) |
| Wash buffer | Same as loading except 10 mM imidazole |
| Elution buffer | Same as loading except 250 mM imidazole |

Ordering Information

| Catalog # | Description |
|-----------------------|---|
| Nickel-Charged | |
| 156-0131 | Profinity IMAC Ni-Charged Resin, 10 ml |
| 156-0133 | Profinity IMAC Ni-Charged Resin, 25 ml |
| 156-0135 | Profinity IMAC Ni-Charged Resin, 100 ml |
| 156-0137 | Profinity IMAC Ni-Charged Resin, 500 ml |
| Uncharged | |
| 156-0121 | Profinity IMAC Resin, 10 ml |
| 156-0123 | Profinity IMAC Resin, 50 ml |
| 156-0125 | Profinity IMAC Resin, 500 ml |
| 156-0127 | Profinity IMAC Resin, 1 L |

Coomassie is a trademark of BASF Aktiengesellschaft.

Profinity IMAC resin is based on UNOsphere bead technology. UNOsphere technology is covered by US patent 6,423,666.

For more information on Profinity IMAC resins, request bulletin 3193.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 2527 3454 Canada (905) 712-2771 China (86 21) 6426 0808 Czech Republic + 420 2 41 43 05 32 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84-0 Greece 30 210 777 4396 Hong Kong (852) 2789 3300 Hungary 36 1 455 8800 India (91-124)-2398112/3/4, 5018111, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 The Netherlands 0318-540666 New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland + 48 22 331 99 99 Portugal 351-21-472-7700 Russia 7 095 721 1404 Singapore 65-64153188 South Africa 00 27 11 4428508 Spain 34 91 590 52 00 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan (886 2) 2578 7189/2578 7241 United Kingdom 020 8328 2000