



Sartobind® IDA 75



The use of a iminodiacetic acid membrane allows purification of proteins via metal ion complex formation.

You may load the metal ion best suited for your application.

Sartobind® IDA provides much higher flow rate than a conventional IDA resin.

Use it with a syringe or connect it via Luer Lock adapter to a liquid chromatography system

Technical Data

Order number	93IDA-42DB-12--V
Quantity	2
Membrane material	Stabilized reinforced cellulose
Pore size	> 3 µm
Number of layers	15
Bed height	4 mm
Bed volume	2.1 ml
Adsorption area	75 cm ²
Ligand	Iminodiacetic acid (IDA)
Coupling method	
Binding capacity for equine heart cytochrome c	> 3 mg/unit
Binding capacity for 6 × His-tagged protein*	7.5 mg/unit
Recommended ions for coupling	Ni ²⁺ , Co ²⁺ , Cu ²⁺ or Zn ²⁺
Flow rate at 0.1 MPa (1 bar, 14.5 psi)	> 10 ml/min
Maximum pressure	0.6 MPa (6 bar, 87 psi)
Housing material	Polysulfone
Chemical stability	Stable in all common chromatography buffers except peroxide and other oxidizing or reactive reagents

* Protected by patents of third parties

**Purification of *E. coli* BL21 (DE3)
pET-bgl-His6-sec***



1 2 3 4 5

Lane 1: Culture supernatant after filtration,
concentrated
Lane 2: Flow-through, concentrated
Lane 3: Wash fraction
Lane 4: Eluate 1
Lane 5: Eluate 2

Loading buffer	0.5 M CuSO ₄
Sample	Supernatant of sonicated <i>E. coli</i> BL21 (DE3) pET-bgl-His6-sec prefiltered with 0.2 µm filter
Washing buffer	50 mM NaH ₂ PO ₄ , 300 mM NaCl, 20 mM imidazole, pH 8.0
Elution buffer	50 mM NaH ₂ PO ₄ , 300 mM NaCl, 250 mM imidazole, pH 8.0
Flow	10 ml/min
Protein eluted	5.9 mg

* Purification of His₆-tagged proteins,
Sartobind Application note SL-4037-e

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA +1.800.3687178
UK +44.1372.737100
France +33.442.845600
Italy +39.055.634041
Spain +34.91.3586102
Japan +81.3.37405407

Sartobind® is a trademark of
Sartorius Stedim Biotech GmbH.

Specifications subject to change
without notice. Printed and copyrighted
by Sartorius Stedim Biotech GmbH
W/sart-199a · G
Publication No.: SL-2041-e07092
Order No.: 85032-533-21