

SARTORIUS

Sartobind® Lab

Think Outside the Column

Incorporating the state-of-the-art in membrane chromatography for affinity capture and charge-based separations, think Sartobind® Lab for rapid antibody capture and purification of most other biomolecules.

Designed to replace legacy resin columns, Sartobind® Lab chromatography units offer:

- Seamless integration into existing processes
- More reliable capture even without a system
- Faster purification in streamlined workflows
- Increased yields and lifetime productivity
- Easier commercialization of biological therapies



Superior Separations

Key Facts

NEW! Sartobind® Rapid A Lab

Rapid capture of antibodies and related molecules for major contaminant reduction.

- <3 min cycles deliver time savings up to 90%
- Resistant to blocking for 100 purifications
- Industry-leading binding capacity ≥35 mg/mL
- Purify 3-17X more antibody per unit

Sartobind® IEX Lab

Faster purification and polishing for most proteins, viruses and nucleic acids.

- <2 min cycles support time savings up to 74%
- Robust matrix capable of 100 purifications
- Dynamic binding capacities up to 29 mg/mL
- Purify up to 7.5X more protein per unit

Ordering Information

| Description | Pack Size | Order No. |
|--------------------------------|-----------|------------------|
| Sartobind® Rapid A Lab, 0.5 mL | 1 | SBLRA025EL-1 |
| | 4 | SBLRA025EL-A |
| Sartobind® D Lab, 2.1 mL | 2 | 93IEXD42DB-12--V |
| Sartobind® Q Lab, 0.4 mL | 4 | 93IEXQ42GB-12--A |
| Sartobind® Q Lab, 2.1 mL | 2 | 93IEXQ42DB-12--V |
| Sartobind® Q Lab, 2.8 mL | 1 | 93IEXQ42BC-12 |
| Sartobind® S Lab, 0.4 mL | 4 | 93IEXS42GB-12--A |
| Sartobind® S Lab, 2.1 mL | 2 | 93IEXS42DB-12--V |
| Sartobind® S Lab, 2.8 mL | 1 | 93IEXS42BC-12 |

Equipment and Accessories

| Description | Pack Size | Order No. |
|-----------------------------------------|-----------|--------------|
| Sartobind® Lab LC system adapter kit | 1 | SBLAAU01-1 |
| Peristaltic pump | 1 | VF-APD0001-1 |
| Peristaltic pump head for 1.6 mm tubing | 1 | VF-APH0001-1 |
| Pump tubing | 1 | VF-ATD0001-1 |

For More Information, Please Contact your Sales Representative