

April 27, 2018

Keywords or phrases:

Redox Cofactors, Plasmids,
Ion Exchange Chromatography,
Membrane Adsorbers, Sartobind® Lab

High Resolution Cofactor and Plasmid Separation by Sartobind® Lab IEX

The phosphate groups on cofactors and the backbone of nucleic acids contribute negative charges to these molecules. They can therefore be easily bound on strong or weak anion exchangers. By use of the included Luer to UNF 10-32 adapters, Sartobind® Lab membrane adsorbers can be connected to most FPLC or HPLC systems for convenient application of complex elution gradients, enabling high resolution separations.

Although the bed height in the three-membrane-layer Sartobind® Lab 15 devices is only 0.8 mm, it is possible to effectively separate separate molecules based on their minimal charge differences, such as the redox cofactors NAD⁺, NADH, NADP⁺ and NADPH, at high resolution (Figure 1).

Sartobind® Lab IEX membrane adsorbers can also be used for the purification of plasmids following alkaline lysis (Figure 2).



Device	Sartobind® Lab Q 15
System	FPLC
Sample	100 µL cofactor mixture (50 µM NAD ⁺ , 50 µM NADH, 150 µM NADP ⁺ , 150 µM NADPH)
Flow Rate	10 mL/min (~25 MV/min)
Buffer A	0.2 mM TEA, pH 7.7
Buffer B	Buffer A + 1 M KCl
Gradient	0% B for 0.8 min 0–20% B in 3.5 min 20–100% B in 1.2 min

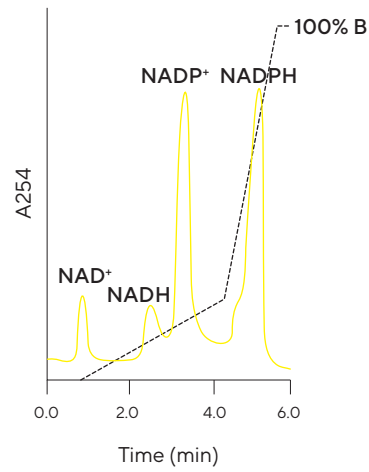


Figure 1: Separation of Cofactors Using Sartobind® Lab Q 15

Device	Sartobind® Lab Q 15
System	FPLC
Sample	1 mL of alkaline lysate containing pBR322, prepared from 100 mL of HB101 culture according to Maniatis et al. (1982)
Flow Rate	0.3 mL/min (~0.75 MV/min)
Buffer A	20 mM Tris pH 8, 1 mM EDTA, 750 mM NaCl
Buffer B	20 mM Tris pH 8, 1 mM EDTA, 850 mM NaCl
Gradient	0% B for 10 min 0–100% B in 20 min

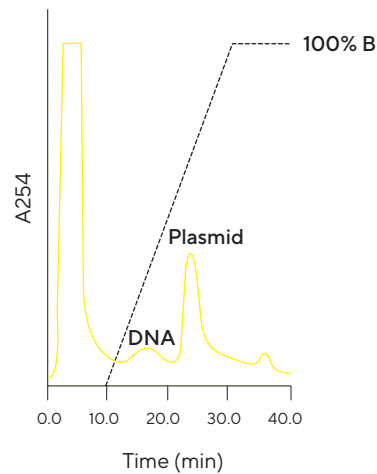


Figure 2: Plasmid Purification Following Alkaline Lysis Using Sartobind® Lab Q 15

Reference

Maniatis, T., Fritsch, E.F. and Sambrook, J. In *Molecular Cloning: A Laboratory Manual* (1982), pp 150–162, Cold Spring Harbor, New York.

Note


Literature published up to c.2022 may reference the use of Sartobind® MA, which is a name previously used for the Sartobind® Lab membrane adsorbers. When these devices were renamed, there was no change made to fit, form or function. Therefore, results collected and methods established using Sartobind® MA devices remain valid also for Sartobind® Lab.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com