

Field Rotor Inspection Program

Protect Your Investment



**BECKMAN
COULTER**

Life Sciences

Save time and get better results

Beckman Coulter offers the widest range of rotors and accessories designed, manufactured and tested as a part of total centrifuge systems. Get the best separation results in the shortest amount of time by using the right rotors. Through 60 years of experience in centrifugation, we have perfected the science of giving you the best equipment for your research needs.



What is the Field Rotor Inspection Program?

The Field Rotor Inspection Program (FRIP) is designed to assist you in maximizing laboratory safety and protecting your investment. FRIP inspections ensure the long life of your laboratory instruments by preventing premature rotor failure due to:

- Stress corrosion
- Metal fatigue
- Wear or damage to anodized coatings

A certified Beckman Coulter Service Specialist will visit your lab and examine all of your rotors.

This includes rotor scoping of all cavities for signs of wear and deterioration. Our specialists will also provide recommendations for replacement of over-speed disks and lubrication of all O-rings and threads as necessary.



A formal training session is also available so that you can learn about proper rotor care and handling, avoiding rotor damage, and the significance of safe and proper laboratory practice.

Centrifuge Rotor Life

The primary factors governing the safe, useful life of any rotor include:

- Expert design and manufacture
- Proper care and handling during use

Expert Design and Manufacture

During the design of a rotor, many tools are used to assure every rotor delivers its expected useful life, is safe, and 100% compatible with approved Beckman Coulter centrifuge systems. Deep expertise in rotor design, computer modeling and simulations, ensures that our rotors support the highest standards. Tests involve many weeks of repeated cycling at maximum speed, and simulations of lab use. These tests often stress the rotors beyond their design speed and expected use. This extensive testing takes safety to the highest level.



Proper Care and Handling

All of our rotors are designed to carry a maximum load at maximum speed. It is important to follow the operational specifications published in each rotor manual since even one run made under conditions of excessive stress may significantly reduce the life of a rotor. Our FRIP experts can help you with care and handling instructions.

Rotor Safety

Each type of rotor has a different useful life based on variables including design life, life-cycle testing and field data. Rotor repair or replacement may be recommended based on these factors, as well as historical use, care and maintenance information. The Field Rotor Inspection Program protects your investment by providing periodic inspections to assess rotor safety and track repair data—ensuring the full, useful lifetime, and best performance of your rotors.



**BECKMAN
COULTER**

***Life* Sciences**

To request your free rotor inspection and learn more about our Field Rotor Inspection Program, visit www.beckmancoulter.com/FRIP

For centrifugation products ordering information, please visit us at:

www.beckmancoulter.com/centrifuge

For Beckman Coulter's worldwide office locations and phone numbers, please visit www.beckmancoulter.com/contact