

TECHNICAL SERVICE & SUPPORT

The “one-stop-shop”
for all things nuclear!



TECHNICAL SERVICE & SUPPORT FROM THE EXPERTS

We are the experts when it comes to difficult process measurements and our technical support team is well known for their level of technical competence, regulatory expertise and commitment to customer service.

All of our service engineers are certified radiation workers and are factory trained.

Please contact us:



+1 (865) 483 1488



berthold-us@berthold.com



Service & Repair Center

Our repair center enables customers to send equipment to Berthold USA to evaluate condition of equipment and determine if it requires and/or is able to be repaired.

Leak Tests / Wipe Tests

Most radioactive materials licenses require sealed sources be leak tested every six months or every three years depending on device registration. We provide do-it-yourself test kits or on-site leak test services.

Nuclear Gauge Shutter Checks

Regulations require operational shutter checks on all devices with shutters, every six months. We can perform this maintenance item for you and identify potential “sticky shutters” before they become a problem.

Nuclear Gauge Commissioning

Commissioning and startup of Berthold devices for General & Specific Licensees.

Nuclear Product Training

Our knowledgeable and experienced engineers can train your personnel to confidently and safely operate, maintain and troubleshoot your Berthold devices.

Preventative Maintenance Plans

Service agreements are available to provide bi-annual leak tests, perform radiation surveys & shutter checks, and audits with training. The scope of work can be tailored to perfectly meet your needs.

Supervisory Services for Installation, Removal and Relocation

Regulatory and radiological safety oversight for plant personnel involved in the physical movement of nuclear gauges.

Source Disposals

Disposal services are not limited to Berthold devices. We can also help you dispose of old Kay Ray, Texas Nuclear, Thermo, Ohmart, Vega, Ronan and other devices. The services include: overseeing the removal of your source from service, providing the required on-site leak testing and radiation surveys, inspecting and packaging the device, preparing for shipment with proper markings, labeling and shipping papers as required by the Dept of Transportation, providing Letter of Transfer of Ownership.



THE EXPERTS IN MEASUREMENT TECHNOLOGY

Berthold Technologies stands for excellent know-how, high quality and superior reliability. The customer is always the focus of our solution. We know our business!

Using our varied product portfolio, our enormous specialized knowledge and extensive experience, we develop suitable solutions together with our customers for new, individual measurement tasks in a wide variety of industries and applications. Berthold Technologies is specialized in radiometric process measurements for 70 years. This is our core competence with state-of-the-art and innovative products and solutions covering a vast range of industries and applications.

We are here for you – worldwide!

The engineers and service technicians from Berthold Technologies are wherever you need them. Our global network assures you fast and above all competent and skilled assistance in case of need. No matter where you are, our highly qualified experts are willing and ready to assist you in a timely manner with the best possible solution for you, even in the most challenging applications..

Berthold Technologies U.S.A., LLC

99 Midway Lane · Oak Ridge · TN 37830
+1 (865) 483 1488 · berthold-us@berthold.com · www.berthold.com/us

