

DETECTOR RETROFIT SOLUTION

Reducing the cost of nuclear source ownership, handling, and disposal



 **BERTHOLD**

DETECTOR RETROFIT SOLUTION TO IMPROVE YOUR PRODUCTION

The challenge

Nuclear measurement sources naturally decay over time, losing their strength, until eventually there isn't enough activity left to reliably measure your process.

Source replacement is costly, time consuming, and does not work towards a reduction in overall site source activity.

In some cases, discontinued products may leave you scrambling to replace failed electronics in a hurry. Fortunately, there is another solution ...

Berthold's solution

Berthold has the most sensitive and stable detectors on the market, a simple replacement allows users to extend the life of their existing sources by years, while improving measurement reliability for all density and level applications.

Berthold detectors can be retrofitted on existing measurements where the source has become too weak to work with the current detector.

With hundreds of systems installed throughout the world, we have the experience and the technical know-how to solve your measurement goals no matter if a new system or a retrofitting solution is required.



Customer benefits

- Extend the life of existing but aged sources
- Managing overall site source activity
- Improving measurement stability and reliability
- Mitigation plan in place well before your measurement quality degrades
- Local support and inventory throughout North and South America

Advantages

- Twice the sensitivity of the competition
- Twice the source life
- Twice the resolution
- Best reliability and stability
- Reduced maintenance costs
- Easy installation on any brand of source holder
- One-time calibration required

High sensitive detectors

Especially in low radiation conditions, our high sensitive detectors improve the signal to noise ratio, what is resulting in superior measurement performance compared to competing systems.





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 · sales-us@berthold-us.com · www.berthold.com

