

ACTIVITY OPTIMIZED SOURCE (AOS)

Optimal mould level control



 **BERTHOLD**

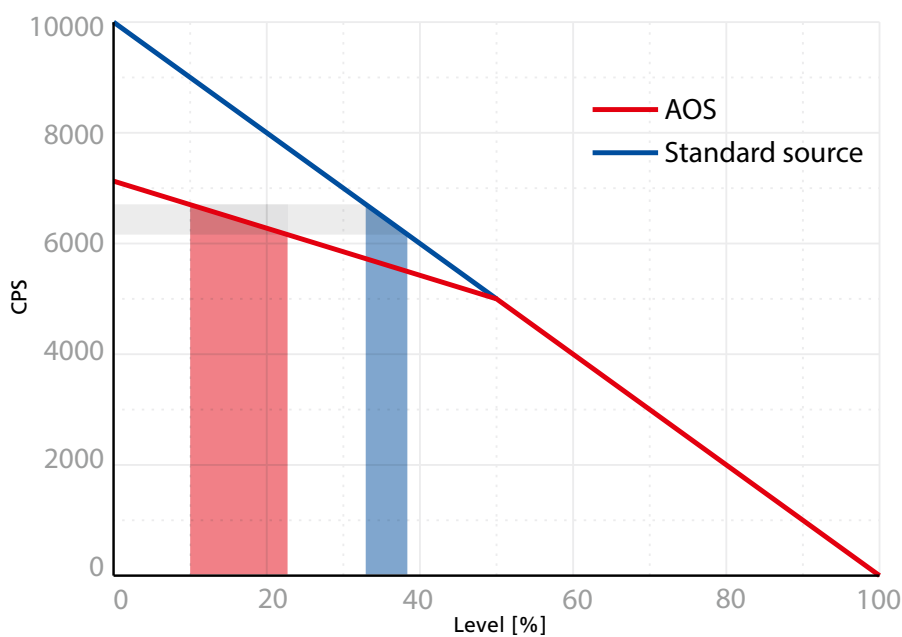
AOS – ACTIVITY OPTIMIZED SOURCE TO IMPROVE YOUR PRODUCTION

The Activity Optimized Source (AOS) is used for the measurement of mould level. The advanced source distribution offers a significant increase of the accuracy in the upper part of the measuring range – where the improved accuracy has very positive effects on the steel quality. The optimized activity distribution accomplishes these enhancements without increasing the total source activity.

Alternatively, when the accuracy is to remain unchanged, the source activity can be reduced by several factors. Especially when installed on small moulds, the AOS is highly efficient in terms of shielding costs and radiation protection matters.

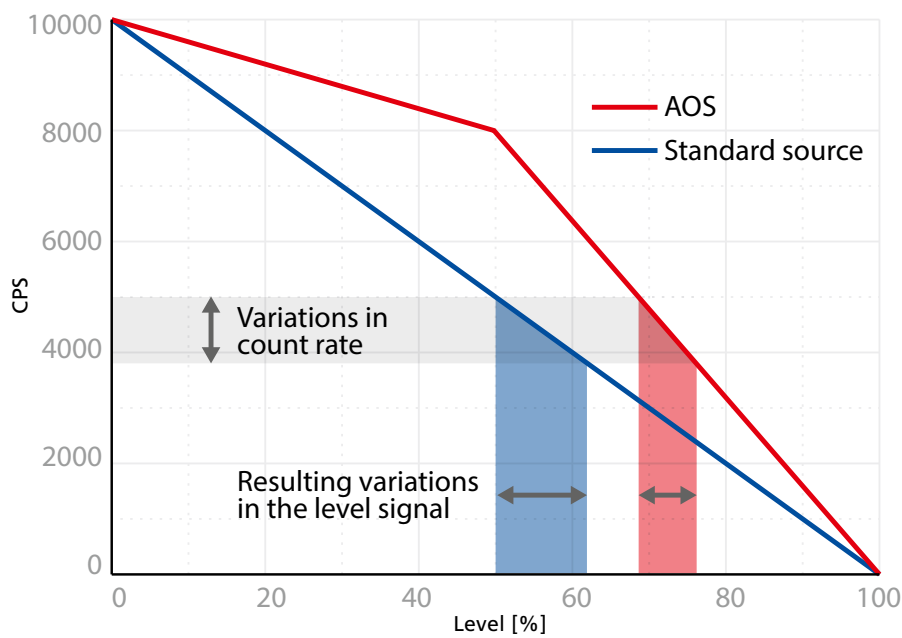
Sources made by Berthold

Berthold operates the only Co-60 rod source production with continuous activity distribution throughout the world. State-of-the-art plant technology is used to produce custom-made sources according to highest quality standards.



Reduced Activity

When using the AOS, the source activity can be significantly reduced, depending on mould size, and measuring conditions. This allows for the use of smaller shielding, which is a major benefit in terms of space and costs. In individual cases it is also possible to work with the cost-effective lead as shielding material instead of Tungsten.



Improved Accuracy

Improved precision at unchanged source activity: This is achieved by changing the activity distribution from the lower part of the source to the upper parts. The lower part of the source is only usable during start of cast and then the accuracy requirement is not so high. The upper part of the source is utilized all the time during normal casting and hence increasing the activity in this region improves casting accuracy. The overall source activity then stays constant.

Features

- Improved precision at unchanged source activity
- Smaller source activity at unchanged precision

Customer benefits

- Best accuracy at the operating point
- Calm level signal short response time
- Small and cost-effective shieldings
- Low dose rates



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 over 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 help needed. 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 GmbH & Co. KG

Calmbacher Straße 22 · 75323 Bad Wildbad · Germany
+49 7081 1770 · industry@berthold.com · www.berthold.com

