

# Confirmation on quality assurance according to nuclear standard KTA 1401

On behalf of the German nuclear power plant operators  
**EnBW Kernkraft GmbH** as a partner of the VGB PowerTech e.V.  
working group "Assessment of Contractors" confirms



## **Berthold Technologies GmbH & Co. KG**

valid for the site

**Calmbacher Straße 22  
75323 Bad Wildbad**



and the scope of supply and services

**Distribution, design, manufacturing and service of  
stationary and handheld radiation measuring  
instruments and systems for radiation protection and  
monitoring in nuclear facilities**



the qualification for system- and product related quality assurance.

The assessment was performed on 10 June 2025 by the firm

**EnBW Kernkraft GmbH**

based on the standard **KTA 1401** as well as on the assessment  
documents of the VGB PowerTech e.V. working group "Assessment of  
Contractors" in consideration of product related requirements.



Details of the assessment are given in the report **0093/20250428/  
0902**.

This confirmation is valid until **01 July 2028** provided that the conditions  
on which the assessment was based have not been changed.



Philippsburg, 10 June 2025

EnKK-RZGDM, Weisenburger

**EnBW Kernkraft GmbH**