

# 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

true for the sites

**Calmbacher Straße 22, 75323 Bad Wildbad  
Gewerbering 35, 76706 Dettenheim**

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 June 14<sup>th</sup>, 2022 by order of EnBW  
Kernkraft GmbH by

**Framatome GmbH**



based on the standard **KTA 1401** as well as on the assessment docu-  
ments of the VGB PowerTech e.V. working group "Assessment of Con-  
tractors" in consideration of product related requirements.



Details of the assessment are given in the report IBGSI/2022/de/0072.

This confirmation is valid until **July 1<sup>st</sup>, 2025** provided that the conditions  
on which the assessment was based have not been changed.

Philippsburg, 2022-06-23

**EnBW Kernkraft GmbH**