

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 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 June 26th, 2019 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/2019/de/0073.

This confirmation is valid until **July 1st, 2022** provided that the conditions
on which the assessment was based have not been changed.



Philippsburg, 2019-07-16

EnBW Kernkraft GmbH