

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

VGB
POWERTECH

Berthold Technologies GmbH & Co.KG Calmbacherstr. 22, D-75323 Bad Wildbad

true for the sites

EnBW

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**

e-on
Kernkraft

the qualification for system- and product related quality assurance.

The assessment was performed on 27th June of 2016 by

EnBW Kernkraft GmbH

RWE
The energy to lead

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 KKP1/2-161/06/16.

Vattenfall Europe
Nuclear Energy

VATTENFALL

This confirmation is valid until **01st of July 2019** provided that the conditions on which the assessment was based have not been changed.

Philippsburg, 28th of June 2016



EnBW Kernkraft GmbH

