



**Nemko-CCL, Inc.**

# Certificate of Compliance

**Certificate:** NA201610530

**Date Issued:** January 20, 2016

**Project:** 257087-7.1

**Issued to:** Berthold Technologies GmbH & Co. KG  
Calmbacher Straße 22  
75323 Bad Wildbad  
Germany

*The products listed below have been certified as being compliant with all applicable requirements of the specifications listed and are eligible to bear the following certification mark*



**Issued by:**

Robert Keller, Senior Engineer/Safety Supervisor

**Authorized by:**

Thomas Jackson, Certification Manager

## **PRODUCTS**

MEASUREMENT, CONTROL, OR LABORATORY EQUIPMENT – Certified to US and Canada Standards

**Product:** Process measurement unit

**Model:** Wall-mounted LB 47x, 1M/3S; Wall-mounted LB 47x, 2M (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety).

**Ratings:** Wall-mounted LB 47x, 1M/3S: 40VA 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x, 2M: 44VA 100-240V, 50/60Hz, Class I



## **APPLICABLE REQUIREMENTS**

UL Std. No. 61010-1 2nd Edition - Safety Requirements for Electrical Equipment for  
Measurement, Control, and Laboratory Use – Part 1: General  
Requirements

CAN/CSA-C22.2 No. 61010-1-04 Second Edition - Safety Requirements for Electrical Equipment for  
Measurement, Control, and Laboratory Use – Part 1: General  
Requirements

This certificate is issued on condition that the holder complies and will continue to comply with the requirements of the above mentioned specifications and  
pursuant to the terms and conditions specified in the Certification Agreement.

## Supplement to Certificate of Compliance

Certificate: NA201610530

Project: 257087-7.1

*Nemko-CCL grants a license to the applicant to apply the Certification Mark to the certified products and that the mark shall only be affixed at the following factory locations*

### Factory Information

Factory Name	Location
Berthold Technologies GmbH & Co. KG	Calmbacher Straße 22 75323 Bad Wildbad Germany

*The products listed, including the latest revision described below, are eligible to be marked in accordance with the referenced Certificate.*

### Product Certification History

Project	Date	Description
257087-7.1	January 20, 2016	<b>Original Certification:</b> <b>Model:</b> Wall-mounted LB 47x, 1M/3S; Wall-mounted LB 47x, 2M (x can be 0 to 8 and describes different software versions for the master and slave modules not affecting safety). <b>Ratings:</b> Wall-mounted LB 47x, 1M/3S: 40VA 100-240V, 50/60Hz, Class I; Wall-mounted LB 47x, 2M: 44VA 100-240V, 50/60Hz, Class I

This Supplement forms an integral part of the Certificate of Compliance

*The certification system, as described in ISO/IEC Guide 67 (Conformity Assessment – Fundamentals of Product Certification), most closely resembles System 3*

Nemko-CCL, Inc. 1940 West Alexander Street Salt Lake City, Utah 84119-2039 Tel (801) 972-6146 Fax (801) 972-8432

