

ISSUE DATE: April 28, 2025

CERTIFICATE AUTHORIZATION NUMBER: 4066



THIS IS TO CERTIFY THAT

Berthold Technologies USA LLC
99 Midway Lane, Oak Ridge, TN 37830

is hereby authorized to continue to apply the
3-A Symbol to the models of equipment, conforming to 3-A Sanitary Standards for:

Number 46-04
46-04 (Refractometers and Energy-Absorbing Optical Sensors)

set forth below

CIP Model: LB 5660-042

VALID THROUGH: **December 31, 2025**

The issuance of this authorization for the use of the 3-A Symbol is based upon the voluntary certification, by the applicant for it, that the equipment listed above complies fully with the 3-A Sanitary Standard(s) designated. Legal responsibility for compliance is solely that of the holder of this Certificate of Authorization, and 3-A Sanitary Standards, Inc. does not warrant that the holder of an authorization at all times complies with the provisions of the said 3-A Sanitary Standards. This in no way affects the responsibility of 3-A Sanitary Standards, Inc. to take appropriate action in such cases in which evidence of nonconformance has been established.

NEXT TPV INSPECTION/REPORT DUE: **April 2030**