



Certificate of Registration

Biotium, Inc.

46117 Landing Parkway, Fremont, CA 94538, USA

have been assessed and approved to:

ISO 9001:2015

Quality Management System

The approved management system applies to the following scope:

Biotium, Inc. located in Fremont, California, manufactures and supplies innovative fluorescence detection technologies and products for life science research, globally. The Quality Management System is applicable to the whole of Biotium, Inc.'s Operations.

Original Approval	21 st October 2014
Current Certificate	18 th February 2023
Certificate Expiry	18 th February 2024
Certificate Number	US3564

Signed: Certification Officer

This certificate remains valid while the holder maintains their administration systems in accordance with the standards and guidelines stated above, which will be audited annually by QAS International. The holder is entitled to display the above registration mark for the duration of this certificate, which should be returned to QAS International upon reasonable request.

Issuing Office: QAS International, 5 Technology Park, Colindeep Lane, London, NW9 6BX, UK

Tel: 020 31989788 • info@qas-international.com • <https://www.qas-international.com>

