



REGISTRATION CERTIFICATE

This document certifies that the administration systems of

Biotium, Inc.

46117 Landing Parkway, Fremont, CA 94538, USA

have been assessed and approved by QAS International to the following management systems, standards and guidelines:

ISO 9001:2015

The approved administration systems apply to the following:

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	21st October 2014
Current Certificate	18th February 2021
Certificate Expiry	18th February 2022
Certificate Number	US3564

Signed: Certification Officer

On behalf of QAS International

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

