



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:

The manufacture and supply of innovative fluorescence detection technologies and products for life science research, globally.

Original Approval:	21 st October 2014
Current Certificate:	18 th February 2025
Certificate Expiry:	18 th February 2026
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

