

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: 21st October 2014

Current Certificate: 18th February 2024

Certificate Expiry: 18th February 2025

Certificate Number: US3564

Signed: Certification Officer

MByas

