



SMITHERS
QUALITY ASSESSMENTS

CERTIFICATE OF APPROVAL

This is to Certify that the Quality Management System of:

AnaSpec, Inc.
34801 Campus Drive
Fremont, CA 94555

has been assessed and approved by Smithers Quality Assessments, Inc., to the following quality management system standards and requirements:

ISO 9001:2015 without Design

The Quality Management System is Applicable to:

Manufacturer/re-packager and supplier of peptides, antibodies, assay kits, synthesis resins, amino acids, dyes, reagents for life sciences research and a distributor of qPCR reagents.

Approval
Certificate Number: 09.357.1

Original Approval: December 23, 2009

Current Certificate: December 12, 2018

Certificate Expires: December 11, 2021



The use of the accreditation mark indicates accreditation in respect of those activities covered by the above certificate number.


on behalf of SQA - J. Michael Hochschwender, CEO

The approval is subject to the company maintaining its system to the required standards which will be monitored by Smithers Quality Assessments, Inc., 121 S. Main St. Suite 300, Akron, Ohio 44308, USA

