

Smithers Quality Assessments, Inc.

CERTIFICATE OF APPROVAL

This is to Certify that the Quality Management System of:

AnaSpec, Inc.
34801 Campus Drive
Fremont, California 94545

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

ISO 9001:2008 without Design

The Quality Management System is Applicable to:

*Manufacturer and supplier of peptides, antibodies, synthesis
resins, amino acids and reagents for life science research.*

Approval
Certificate Number: 09.357.1

Original Approval: December 23, 2009

Current Certificate: December 23, 2009

Certificate Expires: December 22, 2012




on behalf of SQA - J. Michael Hochschwender, President

The use of the accreditation mark indicates accreditation in respect of those activities covered by The Accreditation Certification Number 09.357.1.

The approval is subject to the company maintaining its system to the required standards which will be monitored by Smithers Quality Assessments, Inc., 425 W. Market St., Akron, Ohio 44303-2099

