

STATE OF CALIFORNIA

DEPARTMENT OF PUBLIC HEALTH FOOD AND DRUG BRANCH

DRUG MANUFACTURING LICENSE

AnaSpec, Inc.
2149 O'Toole Avenue
San Jose, CA 95131

LICENSE NUMBER: 41747
EXPIRATION DATE: 11/4/2010

The person named herein is licensed to manufacture drugs through the expiration date of this license. This annual license is issued in accordance with the provisions of Division 104, Chapter 6, Article 6 of the California Health and Safety Code and is not transferable to any other person or place. The licensee is required by law to immediately notify the California Department of Public Health of any change in the information reported in the application.