



| | |
|---|--|
| Revision Number: 1.2 | Last updated December 2013 |
| <u>Product and Company Identification</u> | |
| Product Name: | Anti-Cyclin-E1 (CT), Z-Fish™ |
| Catalog Number: | 55622s |
| Manufacturer/Supplier: | AnaSpec, Inc. 34801 Campus Drive Fremont, CA 94555 Toll-Free: 800-452-5530 Tel: 510-791-9560 Fax: 510-791-9572 http://www.anaspec.com e-Mail: service@anaspec.com |
| <u>Regulation</u> | |
| <p>The "Registry of Toxic Chemical substances" has deemed this material as nonhazardous. Therefore, <u>no material data sheet is required.</u></p> <p>It should be noted that the chemical, physical and toxicological properties of this product have not been investigated thoroughly. Individuals should exercise due care when handling this material.</p> <p><u>This action is in compliance with State and Federal OSHA standards and regulations</u></p> | |
| <u>Other Information</u> | |
| <p><i>This material is sold for research purposes only and is not required to appear on the TSCA inventory. It is not intended for food, drug, household, agricultural or cosmetic use. Its use must be supervised by a technically qualified individual experienced in handling potentially hazardous chemicals. The above information is believed to be correct but does not purport to be all inclusive and shall be used only as a guide. Users should make independent decisions regarding completeness of the information based on all sources available. Anaspec shall not be held liable for any damage resulting from handling or from contact with the above product.</i></p> | |