

## Company Profile



### About AnaSpec

 [www.anaspec.com](http://www.anaspec.com)

#### Integrated Proteomics Solutions

Established in 1993, AnaSpec, Inc. is a leading provider of integrated proteomic and genomic solutions™ for worldwide life science research. As a subsidiary of Kaneka Eurogentec S.A., AnaSpec offers expertise in peptides, detection reagents, antibodies and assay kits. AnaSpec's broad product line of catalog and custom biochemicals and reagents, are used by scientists for basic research, high-throughput screening, and drug discovery.

More than simply providing products or services, AnaSpec is singularly focused on providing solutions for the most demanding research challenges.

AnaSpec is staffed by experts in the fields of peptide technology, fluorescence technology, antibodies, and combinatorial chemistry. The expertise in these complementary fields translates into both premium quality products and cutting-edge innovation.

AnaSpec is certified to ISO 9001:2008. As a company founded by researchers, we recognize the inextricable connection between the quality of one's tools and the success of one's research. Since our inception, AnaSpec has had a longstanding reputation for the unsurpassed quality of our products.

AnaSpec provides unparalleled customer support always ready to answer your order and technical inquiries with personalized, attentive care.

The synergy of AnaSpec's technologies combined with a premium placed on the values of quality, innovation, and service, make AnaSpec uniquely qualified as the leading provider of Integrated Proteomics Solutions™.

### About Kaneka Eurogentec S.A.

 [www.eurogentec.com](http://www.eurogentec.com)

Kaneka Eurogentec is an international company founded in 1985 and part of Kaneka Corp. since 2010. Headquartered in Liege (Belgium) it provides products and services to scientists involved in the life science research, molecular diagnostics and therapeutic developments. Kaneka Eurogentec has production sites in Europe and United States of America. The company is recognized as one of the major suppliers in the field of genomics and proteomics as well as a trusted FDA inspected Contract Development and Manufacturing Organization (CDMO) for the bio-production of pharmaceuticals (vaccines and medicines).

### About Kaneka Corporation

 [www.kaneka.com](http://www.kaneka.com)

Kaneka is an innovation-oriented chemical company. Traditionally the company has been active in polymers, fermentation, biotechnology and electronics, as well as other fields. Business activities now span a broad spectrum of markets ranging from plastics, EPS resins, chemicals and foodstuffs to pharmaceuticals, medical devices, electrical and electronic materials and synthetic fibers. The company has been a pioneer among Japanese chemical companies in establishing overseas operations, beginning in 1970 with a subsidiary in Belgium.