

Instrumentation

All food contaminant AlphaLISA kits were developed on the EnVision multimode plate reader. In addition to fast, sensitive Alpha technology detection, the EnVision Multilabel microplate reader provides fluorescence intensity, luminescence, absorbance, fluorescence polarization, and time-resolved fluorescence detection technologies. The system is based on hybrid technology, combining filters and a monochromator for enhanced flexibility. It incorporates proprietary Direct Double Optics™ technology for high speed and sensitivity in simultaneous filter-based readouts, such as FRET assays, and for TRF, lamp-based excitation is standard, with the option of a high energy laser for higher speed and sensitivity.



Figure 6. EnVision plate reader.

Products and Part Numbers

Table 2. Products for AlphaLISA food contamination detection assays.

AlphaLISA Kit	Size	Part Number
Chloramphenicol AlphaLISA Detection Kit	500 assay points	AL385C
	5,000 assay points	AL385F
Chloramphenicol AlphaLISA Biotin-free Detection Kit	500 assay points	AL393C
	5,000 assay points	AL393F
Aflatoxin AlphaLISA Detection Kit	500 assay points	AL390C
	5,000 assay points	AL390F
Soybean agglutinin AlphaLISA Detection Kit	500 assay points	AL404C
	5,000 assay points	AL404F
Beef troponin AlphaLISA Detection Kit	500 assay points	AL405C
	5,000 assay points	AL405F
½ AreaPlates (96-well, white)	50 plates	6005560
	200 plates	6005569
OptiPlate® (96-well, white)	50 plates	6005290
	200 plates	6005299
OptiPlate (384-well, white)	50 plates	6007290
	200 plates	6007299
AlphaPlate® (384-well, gray)	50 plates	6005350
	200 plates	6005359
AlphaPlate (1536-well, gray)	50 plates	6004350
TopSeal® -A	6050185	