

50+ years of expertise. Applied.

Applied Biosystems™ animal health diagnostics solutions are built on a foundation of proven performance and innovation for the future. They combine decades of research investment and industry expertise—global and local—into a robust, cutting-edge offering from the world leader in serving science.

1949 **Bovine tuberculosis**—production of tuberculin purified protein derivative (PPD) for tuberculosis skin test begins at our facilities in the Netherlands, one of the world's largest PPD production facilities.

1949 **Bovine spongiform encephalopathy (BSE)**—the Prionics™-Check WESTERN kit detects the first BSE-positive cow in routine testing, preventing its meat from entering the food chain and helping avoid human infection.

2004 **Bovine viral diarrhea (BVD)**—the first VetMAX™ real-time PCR kit is developed and launch for the detection of BVD virus in Europe. It's used in European countries to help eradicate the virus and monitor the disease.

2007 **Bluetongue (BT)**—VetMAX™ real-time PCR kits for all BT virus genotypes and kits for subtypes BTV1, BTV2, BTV4, BTV6, BTV8, BTV9, BTV11, and BTV16 detection are approved by the European Community Reference Laboratory (CRL). They are currently used by several European countries for outbreak detection.

2008 **Porcine reproductive and respiratory syndrome (PRRS)**—we launched VetMAX™ PRRSV reagents and controls for RNA detection of NA and EU PRRS virus strains.

2012 **Foot-and-mouth disease (FMD)**—with the addition of the PrioCHECK™ FMDV Type Asia1 and FMDV Type A ELISA test kits to the PrioCHECK™ FMDV NS and Type O kits, we offer the most extensive portfolio of proven solutions for FMD virus detection.

2014 **Avian influenza (AI)**—the VetMAX™-Gold AIV Detection Kit is the first USDA-licensed PCR-based kit for the detection of AI virus. It significantly improves outbreak detection, especially in import/export poultry products.

2014 **Trichomoniasis (Trich)**—the VetMAX™-Gold Trich Detection Kit is the first USDA-approved, real-time PCR test to detect *Trichomonas foetus*. Trich is a sexually transmitted disease that causes embryonic loss and infertility, resulting in significant economic loss within the cattle industry due to open (nonpregnant) and late-calving cows.

2014 **Swine influenza (SI)**—we offer two USDA-licensed VetMAX™-Gold real-time PCR kits for SI virus screening and subtyping to global pork producers for the rapid detection of this highly contagious viral infection of pigs.

2015 **Q fever**—we have the most extensive portfolio of diagnostic PrioCHECK™ ELISA and VetMAX™ real-time PCR kits enabling the detection and quantification of *Coxiella burnetii* in sheep, goats, and cattle. These products also help countries manage Q fever and its zoonotic risk.

Present

Find solutions for current and future challenges at thermofisher.com/animalhealth