



Jena Bioscience, with over 25 years of experience, is a leading provider of innovative and high-quality reagents and customized services in the life science field.



Bulk – Tailored to your scale.

Ready-made or customized life science reagents (liquids, lyo-ready or lyophilized) in large quantities, with personalized packaging and scientific support.



OEM – Your products, your way.

Exclusivity with customized formulations, packaging and branding under strict confidentiality.

Contact our Molecular Biology Experts:

Phone: +49(0)3641 - 6285 000

Mail: molbio@jenabioscience.com



Dr. Bürk Schäfer

Head of Molecular Biology Dept.



Dr. Juliane Buschmann

Business Development Manager



Jena Bioscience

MEAT DETECT qPCR SERIES (HALAL)



Food Control Kits

Jena Bioscience GmbH

Loebstedter Str. 71 – 07749 Jena – Germany

Molecular Biology



MEAT DETECT qPCR SERIES (HALAL)

Pork / Pork+Horse / Pork+Horse+Beef

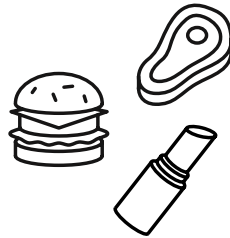
- Most simple handling
- Very fast detection in less than 1 hour
- All required components included

Direct DNA Amplification for Food Control

Sample Collection

Small piece (2-3 mm) of

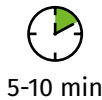
- Raw or cooked meat
- Processed food
- Feed sample
- Cosmetic product



Lysis

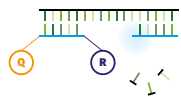


+ Extraction
buffer



5-10 min

Real-time PCR



< 60 min

Cat.-No.

Amount

MeatDetect qPCR Kit Pork (Halal)

PCR-701-24 24 tests

PCR-701-96 96 tests

MeatDetect qPCR Kit Pork+Horse (Halal)

PCR-702-24 24 tests

PCR-702-96 96 tests

MeatDetect qPCR Kit Pork+Horse+Beef (Halal)

PCR-708-24 24 tests

PCR-708-96 96 tests

Pork Genomic DNA

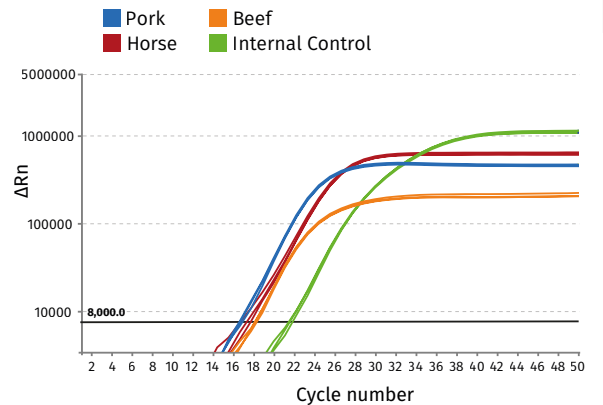
PCR-705 20 µg

Horse Genomic DNA

PCR-706 20 µg

Bovine Genomic DNA

PCR-707 20 µg



MeatDetect qPCR Kit Pork+Horse+Beef

4-plex qPCR amplification plot for a sample with 1 ng DNA each extracted from pork, horse and beef as well as the internal positive control.