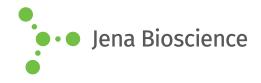
DATA SHEET





Pork Genomic DNA

Positive control template for PCR

Cat. No.	Amount
PCR-705	20 μg

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C, avoid freeze/thaw cycles,

aliquoting of DNA samples is recommended

Shelf Life: 12 months

Form: liquid

Concentration: 200 ng/µl

Description:

Pork genomic DNA is recommended for genomic analysis e.g. as template in positive control PCRs like for the MeatDetect qPCR Kit Pork and MeatDetect qPCR Kit Pork+Horse, #PCR-701 / #PCR-702. The DNA was purified from muscle meat of pork. The DNA fragments have a size of up to 50 kb.

Quality control:

Quality and concentration is confirmed by UV/Vis spectrometry.