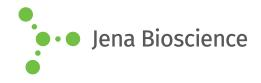
## **DATA SHEET**





## Horse Genomic DNA

Positive control template for PCR

Cat. No.	Amount
PCR-706	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:**

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

## **Quality control:**

Quality and concentration is confirmed by UV/Vis spectrometry.