



## ID-Check Speciation Kits



### Introduction

The identification of meat species present in food and ingredient samples is essential to verify the origin and traceability of raw materials as well as the verification of cleaning processes in production lines. Bio-Rad™ ID-Check Speciation Kits are designed for reliable, specific, and sensitive detection.

ID-Check Speciation Kits are:

- **Available for 8 species** — bovine, chicken, duck, goat, horse, sheep, swine, and turkey meat
- **Ready to use** — user-friendly extraction and detection kits for same-day results in less than 2 hours
- **Reliable** — detects species-specific mitochondrial DNA (mtDNA). High degradation resistance and copy number relative to nuclear DNA (nuDNA) improves detection
- **High sensitivity** — limits of detection at least 0.01%, enabling detection below required limits
- **High specificity** — 100% specificity across multiple species of meat, seafood, fruits, and vegetables, so you can be confident in your results
- **Includes an internal amplification control (IAC)** — improve confidence by excluding false negatives due to PCR inhibition

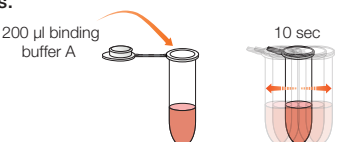
## Easy-to-Use Workflow

ID-Check Speciation Kits are designed for use with Invisorb Spin Tissue Mini DNA Extraction Kits for easy sample preparation and testing.

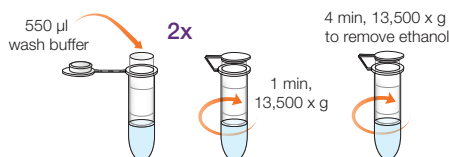
### 1. Lyse samples.



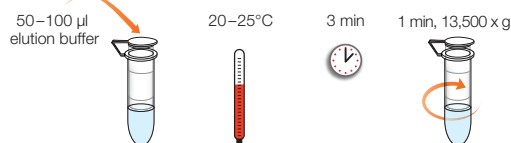
### 2. Bind nucleic acids.



### 3. Wash.



### 4. Elute nucleic acids.



### 5. Run PCR.



## Additional Materials

- Food processor
- Spectrophotometer
- Vortexer
- Microcentrifuge
- Analytical balance
- Thermoshaker (capable of  $52 \pm 1^\circ\text{C}$  and 1,300 rpm)

## Specificity

All ID-Check Speciation Kits have been tested on diverse sample matrices (Table 1) to ensure specificity. No detectable amplification was obtained with any kit for off-target species.

**Table 1. Tested sample matrices for ID-Check Speciation Kits.**

Meat*		
Cow	European wild boar	Rabbit
Chicken	Goat	Sheep
Deer	Horse	Turkey
Duck	Pig	
Fish and Seafood		
Codfish	Grouper	Mussel
Gilt-head bream	Monkfish	Shrimp
Fruits and Vegetables		
Almond	Ginger	Pine nut
Apple	Hazelnut	Pistachio
Apricot	Linseed	Plum
Beans	Lupine	Potato
Brazil nut	Macadamia nut	Pumpkin
Broccoli	Maize	Rice
Cabbage/turnip	Mushroom	Saffron
Carrot	Mustard	Sesame
Cashew	Nutmeg	Soy
Cauliflower	Oats	Sunflower
Celery	Onion	Thyme
Chives	Parsley	Tomato
Cinnamon	Peach	Walnut
Coriander	Peanut	Wheat
Cumin	Pear	White pepper
Garlic	Pecan nut	

\* Depending on the kit being tested.

## Ordering Information

Catalog #	Description
<b>Instruments</b>	
17007991	CFX Opus Deepwell Real-Time PCR System
17007992	CFX Opus 96 Real-Time PCR System
3600037	CFX96 Touch Deep Well Real-Time PCR System
<b>Speciation Kits</b>	
12019385	ID-Check Swine Speciation Kit
12019393	ID-Check Sheep Speciation Kit
12019403	ID-Check Bovine Speciation Kit
12019404	ID-Check Goat Speciation Kit
12019411	ID-Check Duck Speciation Kit
12019421	ID-Check Chicken Speciation Kit
12019423	ID-Check Horse Speciation Kit
12019431	ID-Check Turkey Speciation Kit
<b>Reagent Kit</b>	
12019412	Invisorb Spin Tissue Mini DNA Extraction Kit
<b>Consumables</b>	
17008941	EZ Micro Test Tubes, 1.5 ml graduated micro test tubes
2240110XTU	1.5 ml Conical Tubes, screwcap, sterile

Visit [bio-rad.com/Speciation](https://www.bio-rad.com/Speciation) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Website [bio-rad.com](https://www.bio-rad.com) USA 1 800 424 6723 Australia 61 2 9914 2800 Austria 00 800 00 24 67 23 Belgium 00 800 00 24 67 23 Brazil 4003 0399  
Canada 1 905 364 3435 China 86 21 6169 8500 Czech Republic 00 800 00 24 67 23 Denmark 00 800 00 24 67 23 Finland 00 800 00 24 67 23  
France 00 800 00 24 67 23 Germany 00 800 00 24 67 23 Hong Kong 852 2789 3300 Hungary 00 800 00 24 67 23 India 91 124 4029300 Israel 0 3 9636050  
Italy 00 800 00 24 67 23 Japan 81 3 6361 7000 Korea 82 080 007 7373 Luxembourg 00 800 00 24 67 23 Mexico 52 555 488 7670  
The Netherlands 00 800 00 24 67 23 New Zealand 64 9 415 2280 Norway 00 800 00 24 67 23 Poland 00 800 00 24 67 23 Portugal 00 800 00 24 67 23  
Russian Federation 00 800 00 24 67 23 Singapore 65 6415 3188 South Africa 00 800 00 24 67 23 Spain 00 800 00 24 67 23 Sweden 00 800 00 24 67 23  
Switzerland 00 800 00 24 67 23 Taiwan 886 2 2578 7189 Thailand 66 2 651 8311 United Arab Emirates 36 1 459 6150 United Kingdom 00 800 00 24 67 23

