

Mario Gadanho¹, Sandra Chaves¹, Amanda Manolis², Geoffrey Cottenet³

¹SGS Molecular, Lisboa, Portugal, ²Thermo Fisher Scientific, Microbiology, Basingstoke, UK, ³Nestlé Research Center, Lausanne, Switzerland

1. INTRODUCTION

Food authenticity and fraud are topics of high interest in the food industry and highly controlled by authorities. The complexity of the food supply chain is challenging the abilities of analytical tools used for traceability of ingredients for food production. The most common method to verify species substitution and species identification is Real Time PCR. However, PCR testing is limited by the number of targets that can be simultaneously identified and differentiated. This can be critical, especially when testing highly processed and complex food that often contain multiple different species.

The introduction of Next Generation Sequencing (NGS) into the food sector revolutionizes food authenticity testing. NGS untargeted approach enables accurate detection and differentiation of thousands of different species in each sample using DNA sequencing that is recognized as the most reliable method for species identification.

In the present study an innovative NGS-based approach was tested for meat, fish and spices/herbs species identification.

2. WORKFLOW



- Short DNA fragments suitable for processed food
- 24-36h workflow
- Multiplex and meta-barcoding approach
- In silico analysis for DNA sequence comparison
- Curated Database with thousands of entries

3. RESULTS

3.1 Meat species results

Number of samples tested: 148
Number of species tested: 49
Species mix: up to 3
Spiked level: 1%, 10%, 50%, 100%

<i>Ovis aries</i>	Sheep
<i>Capra hircus</i>	Goat
<i>Lepus capensis</i>	Hare
<i>Oryctolagus cuniculus</i>	Rabbit
<i>Macropus rufus</i>	Kangaroo
<i>Capreolus capreolus</i>	Roe Deer
<i>Cervus elaphus</i>	Red Deer
<i>Rangifer tarandus</i>	Reindeer
<i>Antidorcas marsupialis</i>	Springbok
<i>Equus hemionus</i>	Zebra
<i>Lama glama</i>	Lama
<i>Gallus gallus</i>	Chicken
<i>Canis familiaris</i>	Dog
<i>Bison bison</i>	Bison
<i>Cervus dama</i>	Fallow Deer
<i>Equus caballus</i>	Horse
<i>Sus scrofa</i>	Pork
<i>Bos taurus</i>	Beef
<i>Meleagris galopavo</i>	Turkey
<i>Cairina moscata</i>	Duck
<i>Alopochen aegyptiacus</i>	Goose
<i>Struthio camelus</i>	Ostrich
<i>Columba livia</i>	Pigeon
<i>Numida meleagris</i>	Guinea fowl
<i>Dromaius novaehollandiae</i>	Emu

<i>Tragelaphus strepsiceros</i>	Kudu
<i>Felix catus</i>	Cat
<i>Rattus norvegicus</i>	Rat
<i>Vulpes vulpes</i>	Fox
<i>Alces alces</i>	Elk
<i>Coturnix japonica</i>	King quail
<i>Bubalus bubalis</i>	Buffalo
<i>Camelus dromedarius</i>	Camel
<i>Crocodylus niloticus</i>	Crocodile
<i>Lophura inornata</i>	Pheasant
<i>Oryx leucoryx</i>	Oryx gazella
<i>Alcelaphus buselaphus</i>	Gnu
<i>Bos grunniens</i>	Cattle Yak
<i>Equus asinus</i>	Donkey
<i>Meles meles</i>	Badger
<i>Tragelaphus scriptus</i>	Antelope
<i>Corvus macrorhynchos</i>	Daw
<i>Mustela erminea</i>	Weasel
<i>Ondatra zibethicus</i>	Muskrat
<i>Anas sp</i>	Mallard duck
<i>Crocodylus siamensis</i>	Crocodile
<i>Phasianus colchicus</i>	Pheasant
<i>Alectoris chukar</i>	Partridge
<i>Aepyercos melampus</i>	Impala

3.2 Plants species (spices/cereals) results

Number of samples tested: 347
Number of species tested: 39
Species mix: up to 5
Spiked level: 1%, 5%, 10%, 20%, 100%

<i>Origanum sp.</i>	Origanum
<i>Allium schoenoprasum</i>	Wild chives
<i>Allium sativum</i>	Garlic
<i>Anethum graveolens</i>	Dill
<i>Argemone sp</i>	Prickly poppy
<i>Avena sativa</i>	Oat
<i>Brassica napus</i>	Rape
<i>Capsicum annum</i>	Cayenne pepper
<i>Carum carvi</i>	Caraway
<i>Cerantonia siliqua</i>	St John's bread
<i>Conium maculatum</i>	Poison henlock
<i>Coriandrum sativum</i>	Coriander
<i>Crocus sativus</i>	Saffron
<i>Cuminum cyminum</i>	Cumin
<i>Curcuma longa</i>	Turmeric
<i>Elettaria cardamomum</i>	Cardamom
<i>Foeniculum vulgare</i>	Sweet fennel
<i>Glycine max</i>	Soybean
<i>Hordeum vulgare</i>	Barley
<i>Juniperus communis</i>	Juniper

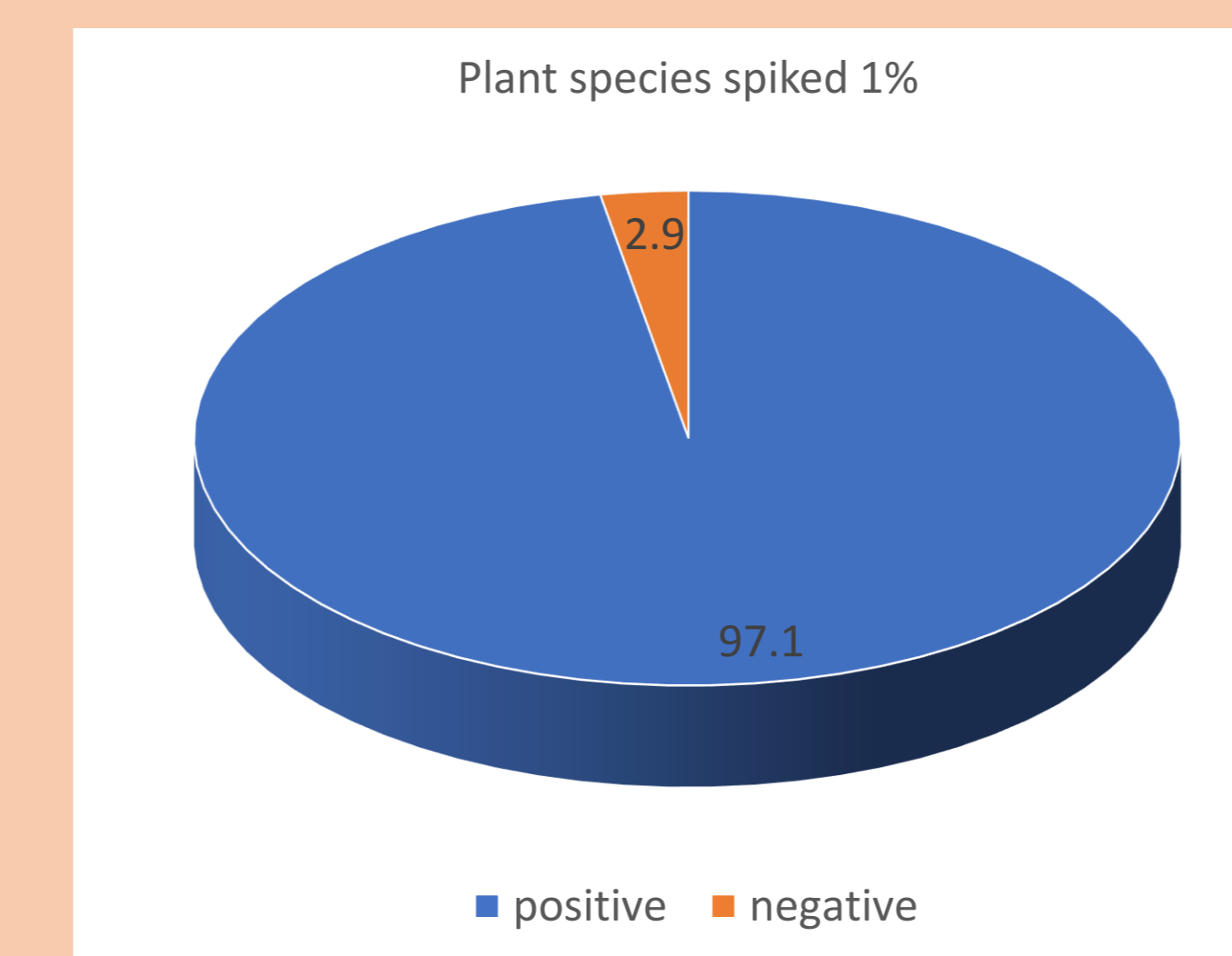
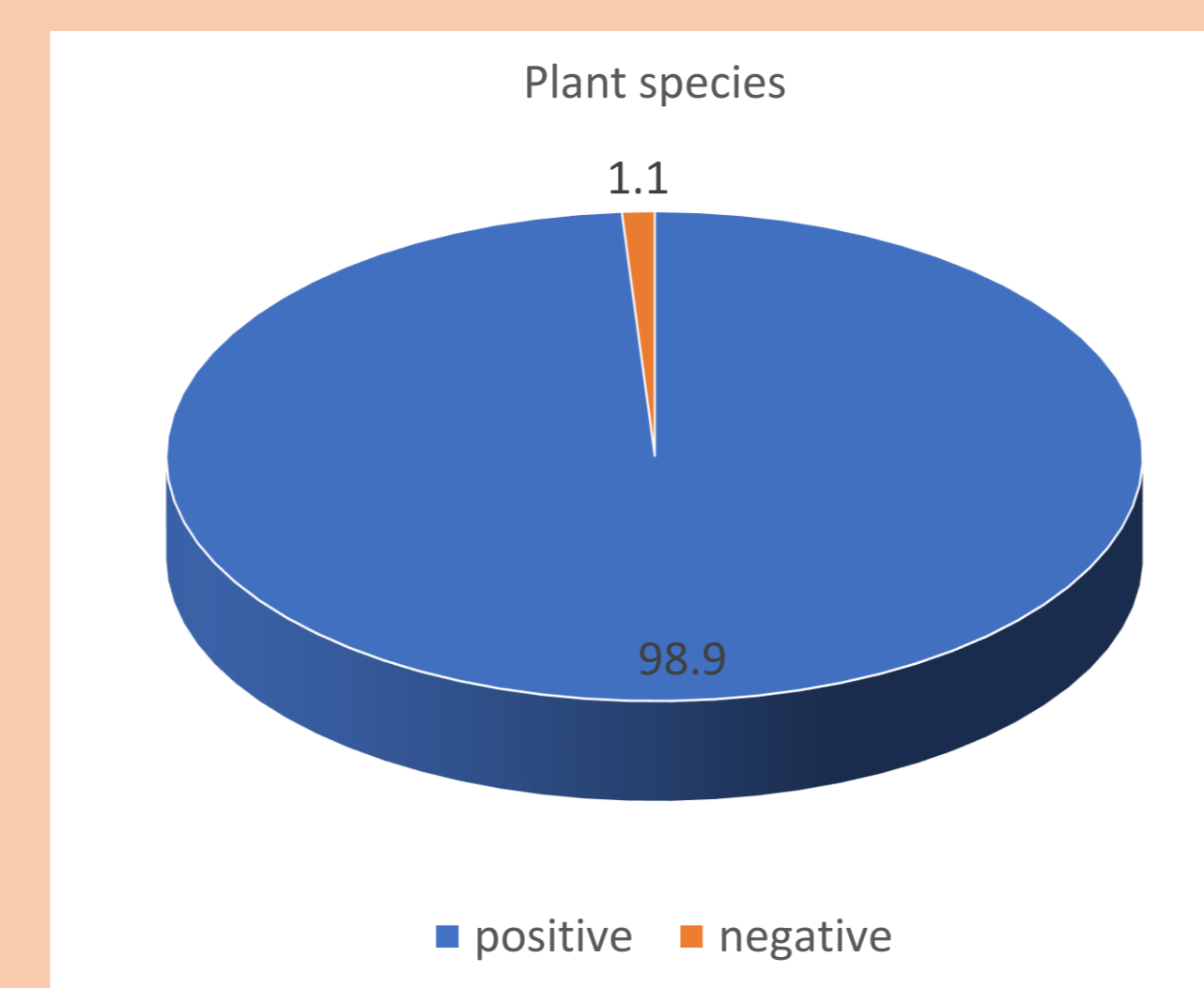
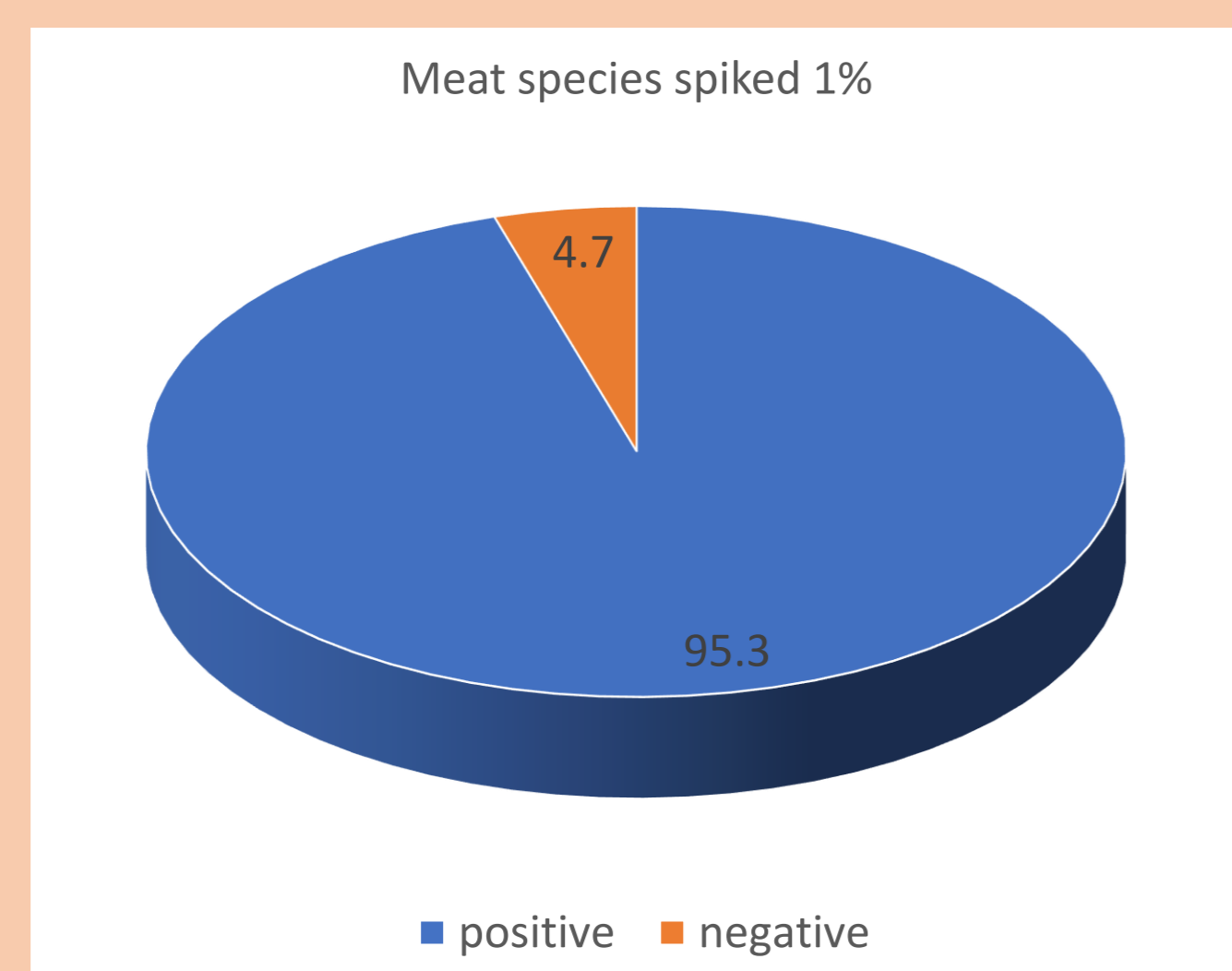
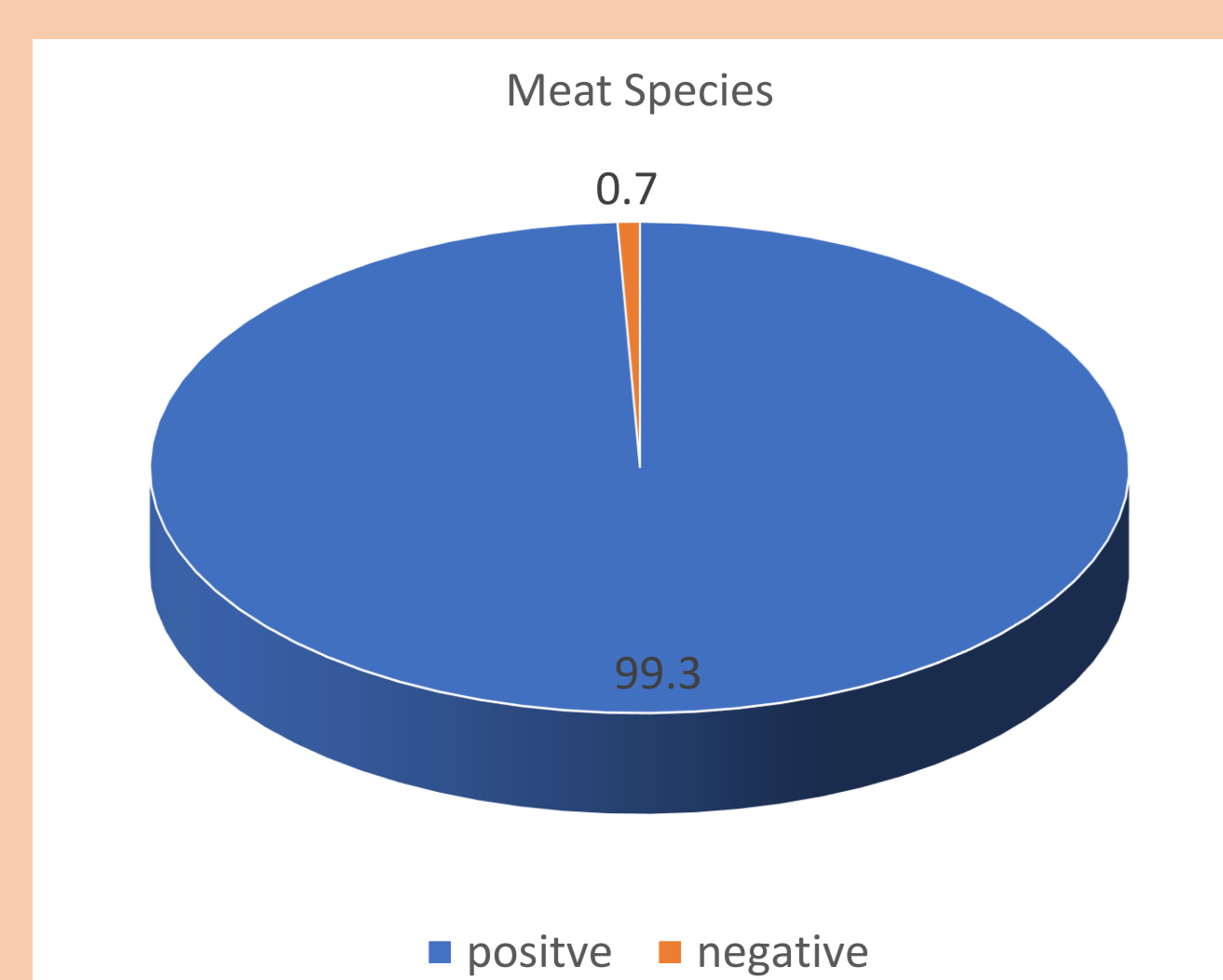
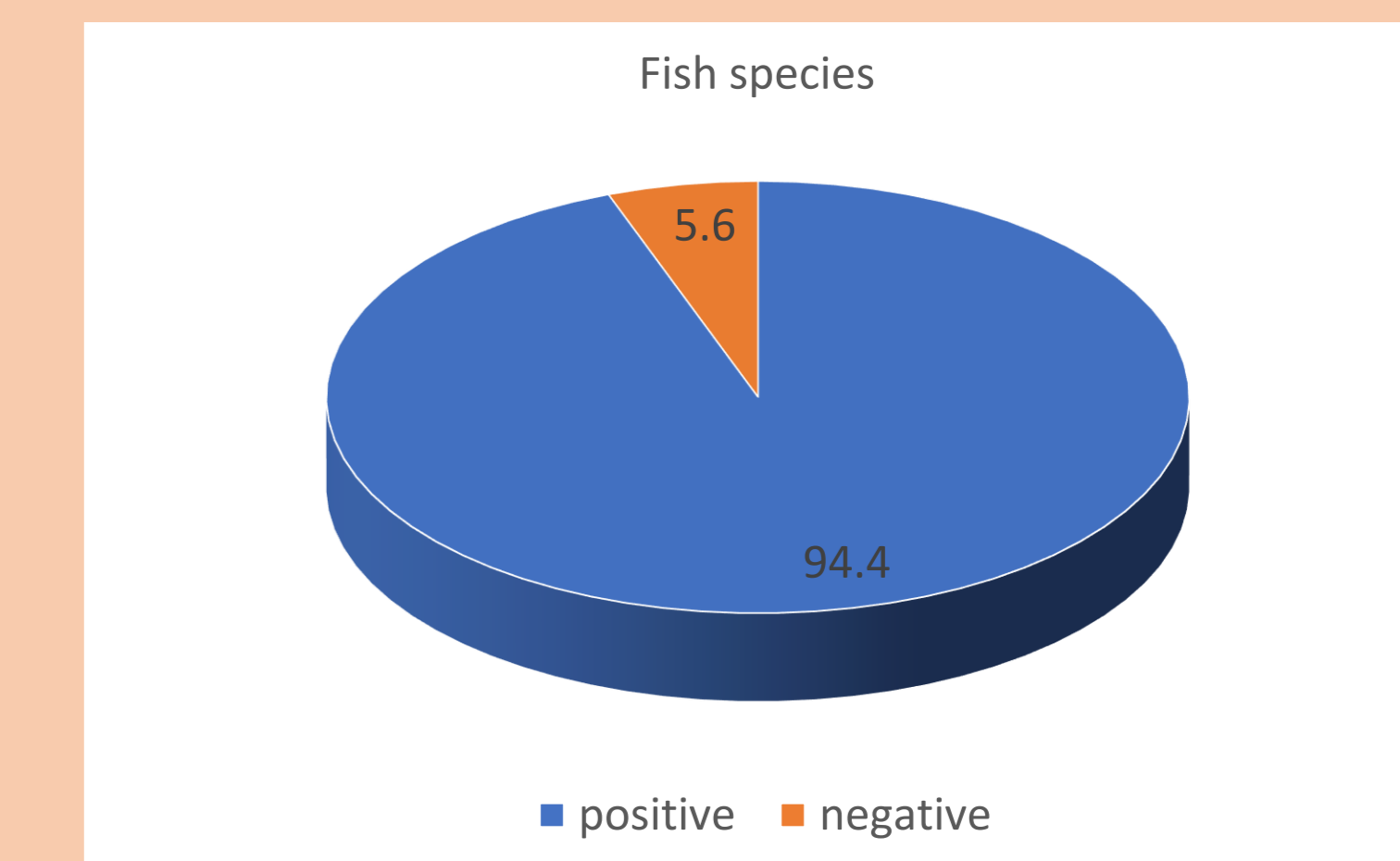
<i>Laurus nobilis</i>	Sweet bay
<i>Manihot esculenta</i>	Cassava
<i>Mentha spicata</i>	Spearmint
<i>Myristica fragrans</i>	Nutmeg
<i>Ocimum basilicum</i>	Sweet basil
<i>Oryza sativa</i>	Rice
<i>Panicum miliaceum</i>	Millet
<i>Papaver somniferum</i>	Opium poppy
<i>Petroselinum crispum</i>	Parsley
<i>Pimpinella anisum</i>	Anis
<i>Piper nigrum</i>	Black pepper
<i>Rosmarinus officinalis</i>	Rosemary
<i>Sesamum indicum</i>	Sesame
<i>Sinapis alba</i>	White mustard
<i>Sorghum bicolor</i>	Sorghum
<i>Thymus vulgaris</i>	Garden Thyme
<i>Triticum aestivum</i>	Wheat
<i>Triticum durum</i>	Durum wheat
<i>Zingiber officinale</i>	Garden Ginger

3.3. Fish species results

Number of samples tested: 78
Number of species tested: 26
Species mix: up to 2
Spiked level: 1%, 2%, 5%, 10%, 100%

<i>Salmo salar</i>	Atlantic Salmon
<i>Thunnus albacares</i>	Yellowfin tuna
<i>Gadus morhua</i>	Atlantic cod
<i>Hippoglossus hippoglossus</i>	Pacific halibut
<i>Limanda limanda</i>	Common dab
<i>Merluccius merluccius</i>	European hake
<i>Melanogrammus aeglefinus</i>	Haddock
<i>Katsuwonus pelamis</i>	Skipjack tuna
<i>Thunnus alalunga</i>	Albacore
<i>Pleuronectes platessa</i>	European plaice
<i>Molva molva</i>	Ling
<i>Sander lucioperca</i>	Pike-perch
<i>Pollachius pollachius</i>	Pollack

<i>Trisopterus luscus</i>	Norway pout
<i>Cynoglossus senegalensis</i>	Witch flounder
<i>Oncorhynchus chrysogaster</i>	Pink salmon
<i>Lophius piscatorius</i>	Angler
<i>Oncorhynchus nerka</i>	Sockeye salmon
<i>Pangasianodon hypophthalmus</i>	Silver carp
<i>Scorpaenopsis diabolus</i>	Atlantic mackerel
<i>Oncorhynchus gorbuscha</i>	Pink salmon
<i>Merluccius hubbsi</i>	Argentine hake
<i>Merluccius productus</i>	North Pacific hake
<i>Macruronus magellanicus</i>	Patagonian grenadier
<i>Merluccius gayi</i>	South Pacific hake
<i>Thunnus obesus</i>	Bigeye tuna



4. CONCLUSIONS

- The NGS workflow could enable to identify all the species tested individually
- Identification of pure material at spiked levels higher than 5% was successfully obtained for more than 94% of the samples tested. A few negative results were obtained for some highly processed food products and mostly for fish-based samples.
- At the level of 1% spiking, more that 95% of the samples were correctly assigned to the species name.
- When combined, all targets could be analyzed simultaneously in a single NGS run which reduce significantly the NGS costs
- Based on these invitro analysis and on the in silico analysis of DNA sequences a database containing thousands of species was created.
- Extension of the databases is an on-going activity to include additional species