

# EXAMPLES of FOOD TYPES ANALYSED

## MEATS

PORK MINCE\*  
 BEEF MINCE\*  
 TURKEY MINCE  
 BEEF SAUSAGE  
 CHICKEN  
 COOKED HAM  
 COOKED CHICKEN  
 PEPPERED SALAMI  
 PRAWNS  
 FISH  
 PORK SAUSAGE  
 SPANISH CHORIZO  
 LASAGNE  
 GERMAN SALAMI  
 MEAT PIES  
 PET FOOD  
 BEEF TRIM

\*Many Retailer "own brands" tried

## DAIRY

MILK  
 YOGURTS  
 SOFT CHEESES  
 HARD CHEESES

## SPICES & NUTS

NUTMEG  
 PEPPER CORNS  
 CASHEW NUTS  
 COCONUTS  
 WALNUTS

## OTHERS

SEAFOOD  
 CHOCOLATE  
 COCO POWDER  
 READY MEALS  
 PIZZA  
 XANTHAM GUM

## PRODUCE

LETTUCE  
 SPINACH  
 VEGETABLES  
 FROZEN VEGETABLES  
 FRESH FRUIT  
 EGG

## JUICES

ORANGE JUICE  
 TOMATO JUICE  
 PINEAPPLE JUICE  
 APPLE JUICE

## ENVIRONMENTAL

COTTON BUD SWABS  
 FECES  
 LITTERS (CHICKEN)