

X-ray detection

# Enhancing performance & efficiency

## NextGuard X-ray inspection new application solutions

Find metallic and non-metallic foreign objects and eliminate “wet” product effect common with metal detectors with the Thermo Scientific™ NextGuard™ X-ray Detection Systems. Designed for a wide variety of applications, NextGuard systems offer enhanced capabilities to inspect packaged products for missing pieces or components, under and over-fills, and other quality problems with easily customizable vision software.

Proper placement of product on the conveyor when passing through the inspection tunnel is imperative in many applications. Some products moving down the production line may need to be “guided” into the center of the conveyor prior to entering x-ray inspection systems. Other small/narrow and/or light weight products can pose a challenge in passing through the shielding curtains into and out of the x-ray tunnel. The following are solutions to solve these challenges for manufacturers.

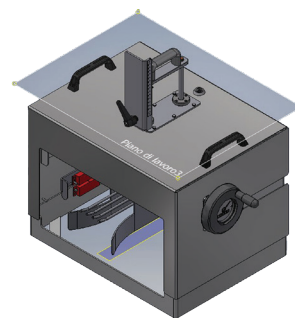


### Adjustable guide rails for single or multi-lane inspection

**Product:** For use on NextGuard C330

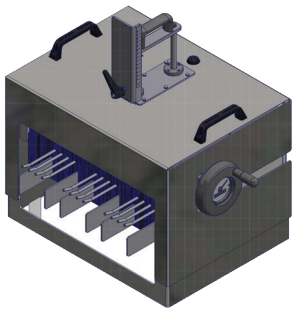
**Application solution:**

- Special radiation shield equipped with adjustable guiding rails – integrated into aperture
- Allows for two-dimensional (vertical & lateral) manual adjustment
  - Run different sized products on the same line



Single lane guide rails

- **Single lane** - helps to guide products into the center area of the inspection tunnel
  - Ensure correct positioning for next step in manufacturing process (i.e. case packing)
  - Aligns product to optimize mass estimation, product verification, or edge masking applications



**Multi-lane guide rails**

- **Multi-lane** - allows for expansion (up to three lanes) of inspection capacity
  - Used in conjunction with multi-lane packaging equipment
  - Guiding products into two or three lanes prior to entering aperture

**Additional applications:** Food & Beverage, Pharmaceutical, Personal Care products



## Thin slit curtains

**Product:** For use on NextGuard C330

**Application solution:**

- Light weight, low grip
- Lead free, FDA approved material – can come into contact with unpackaged food items
- Curtain strips are 4mm wide
- Use with low profile/narrow, light weight products

**Additional applications:** Food & Beverage, Pharmaceutical, Personal Care products



Contact us today [thermofisher.com/productinspection](https://www.thermofisher.com/productinspection)

**thermo**scientific