## thermo scientific

TRUNARC IN ACTION

## Protecting officers in the field

## Porter County Sheriff's Office

Concerned their officers were at risk due to fentanyl exposure, the Porter County Sheriff's Office deployed the Thermo Scientific<sup>™</sup> TruNarc<sup>™</sup> Handheld Narcotics Analyzer. Officers can scan suspected narcotics in the packaging, without having to directly handle them. This significantly lowers officer risk of exposure.

- Previously, officers had to physically remove a sample of the substance from its bag
- Officers could inhale or come in contact with substances that could be extremely dangerous
- The digital library on TruNarc can be updated as new threats emerge
- The device identifies precursors and cutting agents, such as baking soda
- Results reports are generated and can be used for court submission



"We are very fortunate that we have this available for us and our officers. Given what's in drugs today, it's much safer for our officers."

 Sgt. Jamie Erow, Public Information Officer for the Porter County Sheriffs Office

## Find out more at thermofisher.com/trunarc

