thermoscientific

TRUNARC IN ACTION

Testing for over 250 narcotics, including fentanyl in a single test

Charles County Sheriff's Office

When officers discovered narcotics on the scene, identifying them was a time consuming and risky exercise. Officers would often handle the drugs as they attempted to determine what the drug was. A small quantity of fentanyl coming in contact with an officer can result in an overdose or death.

- The sheriff's office received a grant to purchase a Thermo Scientific™ TruNarc™ analyzer
- One scan of the sample can test it against all of the narcotics in the TruNarc library
- Samples can be tested without removing them from the plastic bag, drastically reducing risks of exposure for officers
- By quickly identifying the presence of fentanyl mixed with heroin officers are able to ensure officer safety when handling the situation
- If a suspect has overdosed, by quickly determining what the drug involved is, officers can get the best treatment and result for the individual
- The TruNarc results can be used to press charges against the suspect immediately instead of waiting months for lab results to come back



"The safety of our officers is one of our top priorities, and the TruNarc allow us to safely identify controlled dangerous substances quickly and accurately without having to wait on lab results. The community will benefit because we will be able to charge suspects sooner and more precisely."

- Charles County Sheriff's Office

