

Misread labels lead to sample mix-up

Dropped samples lead to waste



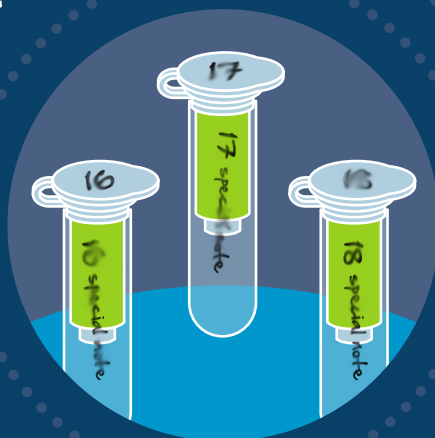
Is that a 6 or a 9?



Numerous pipetting steps require excess focus



Smeared sample labels cause confusion



Back and forth and back and forth



# Alleviate Your Pain Points by Taking EMnetik for a Spin

Nucleic acid cleanup is ripe for innovation. For decades, column cleanup has required excessive time and resources from labs. What's more, the procedures are prone to user error. It's time to bring nucleic acid cleanup into the 21st century.



When paired together, the **EMnetik 24** and the **EMnetik PCR** clean up kit are a one-stop-shop for nucleic acid cleanup. The innovative bead technology and new micro-particle processor streamlines and expedites column cleanup procedures, requiring less time and fewer touchpoints while performing the entire nucleic acid cleanup process in a single space. With EMnetik 24 and its accompanying products, nucleic acid cleanup is no longer a burden. Now, you have time to free your mind.

