

CHANGE THE WAY
YOU WORK
CHANGE THE WAY
WE LIVE

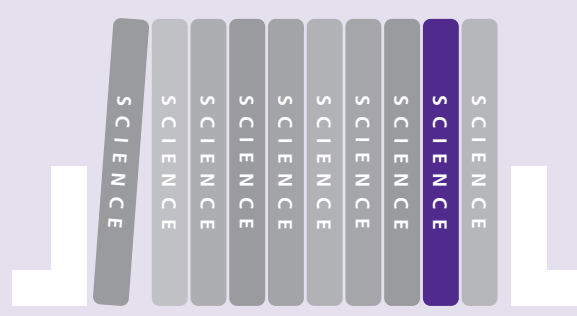
Automate your flame AA sample introduction for unparalleled productivity, reproducibility, and profitability.

THE PROBLEM

14,350 
ILLNESSES FROM
FRESH PRODUCE
BETWEEN
1996 AND 2010

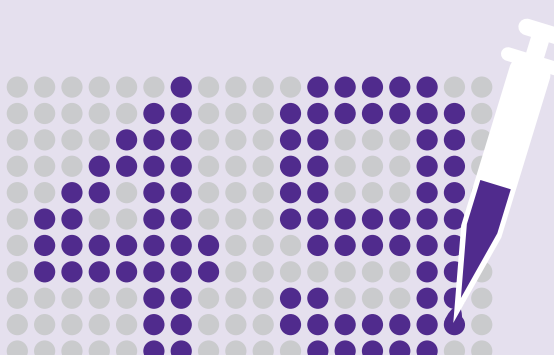

3M Tons of toxic
chemicals discharged
into the air, water
and soil each year

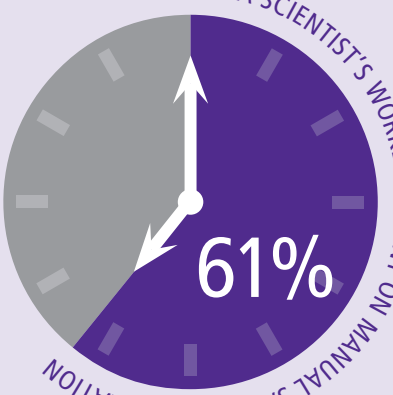
THE CHALLENGES


ONLY **10%** OF PUBLISHED
APPLICATIONS ARE REPRODUCIBLE

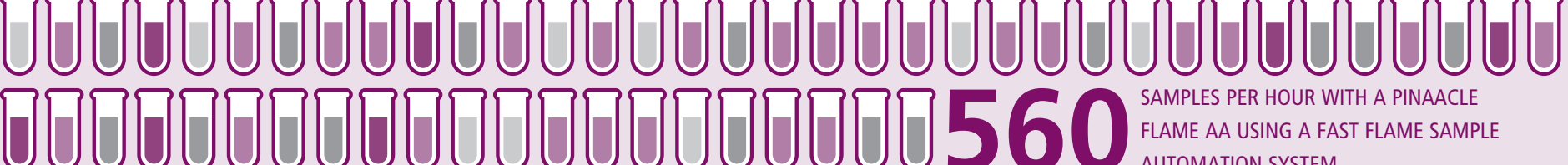

AVERAGE NUMBER OF COUNTRIES SUPPLYING
INGREDIENTS FOR A TYPICAL AMERICAN MEAL:


5




PERCENT OF OUT-OF-SPECIFICATION
ERRORS CAUSED BY
MANUAL SAMPLE PREP


61% OF A SCIENTIST'S WORKDAY SPENT ON MANUAL SAMPLE PREPARATION

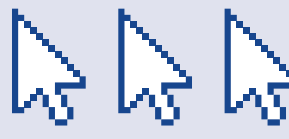
THE SOLUTION


560 SAMPLES PER HOUR WITH A PINACLE
FLAME AA USING A FAST FLAME SAMPLE
AUTOMATION SYSTEM  = 10 SAMPLES

FAST Flame is nearly
3X faster
than other automated systems

 **5mL** = \$\$\$
SAMPLE USED LESS CONSUMPTION
AND REDUCED
OPERATING COSTS 

THE BENEFITS

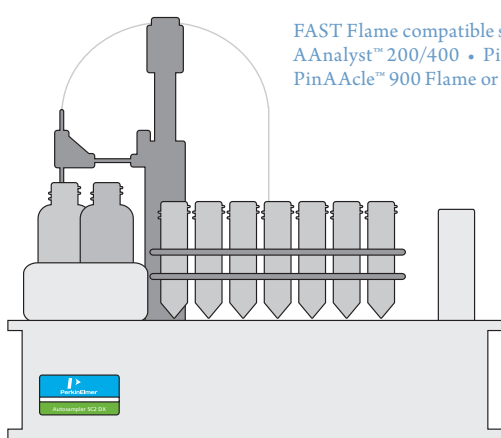

3 CLICKS
FROM SETUP TO ANALYSIS
USING SYNGISTIX™ FOR
AA SOFTWARE


RISK %
NO RISK OF HUMAN ERROR


Syngistix Enhanced Security™ option
simplifies 21 CFR Part 11 compliance


50 µg/L
DETECTION LIMITS

 **2 17 00**
ONE THOUSAND REPLICATE ANALYSES IN TWO HOURS SEVENTEEN MINUTES



FAST Flame compatible systems:
AAAnalyst™ 200/400 • PinAAcle™ 500 Flame AA
PinAAcle™ 900 Flame or Flame/Graphite Furnace AA Systems

FAST Flame Sample Automation Platform for flame AA applications—www.perkinelmer.com/FASTflame



For the Better